



Nonradioactive Iodide Assay kit

A brand name of

bertin
pharma

For laboratory research use only. Not for human or veterinary diagnostic use.

Bertin Pharma also markets pre-analytical products, EIA kits, antibodies & biochemicals for:

- ▶ **Cardiology / Hypertension**
- ▶ **Diabetes / Obesity**
- ▶ **Endocrinology / Metabolism**
- ▶ **Inflammation**
- ▶ **Pharmacology**
- ▶ **Psychopharmacology**
- ▶ **Nitric Oxide**
- ▶ **Oncology / Apoptosis**
- ▶ **Oxidative injury**
- ▶ **Cell signaling**
- ▶ **Drug metabolism**

Do not hesitate to contact our after-sales services for further information at bioreagent@bertinpharma.com



**Nonradioactive Iodide Assay kit
#D05076.96 wells**

For research laboratory use only
Not for human diagnostic use

This assay has been developed & validated
by Bertin Pharma



Fabriqué en France
Made in France

#D11076
Version: 0117

Table of contents

| | |
|--|----|
| ▶ Precaution for use | 6 |
| ▶ Iodide background | 8 |
| ▶ Principle of the assay | 10 |
| ▶ Materials and equipment required | 11 |
| ▶ Sample collection and preparation | 12 |
| ▶ Reagent preparation | 13 |
| ▶ Assay procedure | 16 |
| ▶ Data analysis | 19 |
| ▶ Acceptable range | 20 |
| ▶ Typical results | 21 |
| ▶ Assay validation and characteristics | 22 |
| ▶ Troubleshooting | 23 |
| ▶ Bibliography | 24 |

96 wells
Storage: +4°C
Expiry date: stated on the package

This kit contains:

| Designation | Item # | Quantity per kit | Form |
|---------------------------------------|----------------|------------------|--------|
| Covered 96 well Microtiter plate | D08000.1 ea | 1 | - |
| Nonradioactive iodide assay reagent A | D22076.100 dtn | 1 | Liquid |
| Nonradioactive iodide assay reagent B | D20076.100 dtn | 1 | Liquid |
| Nonradioactive iodide assay standard | D06076.1 ea | 1 | Liquid |
| Technical Booklet | D11076.1 ea | 1 | - |
| Well cover sheet | - | 1 | - |

Each kit contains sufficient reagents for one 96-well plate. This allows for the construction of one standard curve in duplicate and the assay of 37 samples in duplicate.

▶ **Precaution for use**

Users are recommended to carefully read all instructions for use before starting work.

Each time a new pipette tip is used, aspirate a sample or reagent and expel it back into the same vessel. Repeat this operation two or three times before distribution in order to equilibrate the pipette tip.

- > For research laboratory use only
- > Not for human diagnostic use
- > Do not pipet liquids by mouth
- > Do not use kit components beyond the expiration date
- > Do not eat, drink or smoke in area in which kit reagents are handled
- > Do not use organic solvent (DMSO, EtOH, MeOH, MeCN, acetone)
- > Avoid splashing

The total amount of reagents contains less than 23 mg of arsenic(III) oxide.

Wearing gloves, laboratory coat and glasses is recommended when assaying kit materials and samples.



Waste elimination:

These reagents are dangerous for environment.

- Do not throw in the sink.
- Use appropriate recovery can.

▷ **Temperature**

Unless otherwise specified, all the experiments are done at room temperature (RT), that is around +20°C. Working at +25°C or more affects the assay and decreases its efficiency.

▶ Iodide background

Present in nature in the form of iodide and iodate, iodine is a solid halogen at normal temperature. It is used in medicine, in the pharmaceutical and food industry. Food is the principal daily supply of iodide in human body [1].

Iodide is important in basal metabolism and permits temperature regulation, intellectual development for children, muscular development, normal heart function and growth of skeleton. Iodide transport is the basis for an emerging approach of selective cancer cell destruction. [2-3]

Iodide uptake from blood into thyroid follicular cells is the first step in the biosynthesis of thyroid hormones T4 and T3, known to regulate many essential biological processes [4-5].

Thyroid hormones are indispensable for body development. This transport is mediated by NIS (sodium iodide symporter), an intrinsic membrane glycoprotein located in the basolateral membrane of thyrocytes.

Since the discovery of NIS, thorough biochemical analysis has elucidated the mechanism of basolateral iodide transport and revealed the key role of NIS in thyroid diseases such as thyroid cancer, autoimmune disease, and congenital hypothyroidism [6].

If rate is not in the normal proportion, some diseases can be developed as underactive thyroid if the rate is too down or overactive thyroid if the rate is too up. Other diseases exist as

chronic thyroiditis of Hashimoto or cancer of the thyroid gland **[7]**.

Iodide deficiency is at origins of many thyroid metabolism disorders, this is why it is important to control rate of iodide to prevent all of these diseases.

▶ Principle of the assay

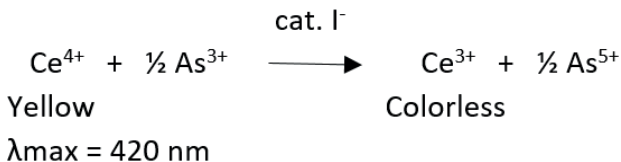
The present assay is a nonradioactive method for the measurement of iodide.

This Iodide Assay is based on the oxido-reduction reaction: cerium(IV) is reduced by arsenic(III). The reduction of yellow (420 nm) cerium(IV) to colorless cerium(III) by arsenic(III) proceeds very slowly but traces of iodide strongly accelerate this reaction with the rate being directly proportional to iodide concentration.

For a given time, decay is inversely proportional to iodide concentration in well.

This method is simple and nonradioactive, and as such it can be used widely.

The principle of the assay is summarized below:



► **Materials and equipment required**

In addition to standard laboratory equipment, the following material is required:

- > Precision micropipettes (20 to 1000 μL)
- > Spectrophotometer plate reader (414 nm or 420 nm filter)
- > Orbital microplate shaker
- > Multichannel pipette and disposable tips 30-300 μL
- > UltraPure water #A07001.1L
- > Polypropylene tubes



Water used to prepare all diluted reagents must be UltraPure (deionized & free from organic contaminant traces).

Otherwise, organic contamination can significantly affect the assay.

Do not use distilled water, HPLC-grade water or sterile water.

UltraPure water may be purchased from Bertin Pharma: item #A07001.1L.

▶ **Sample collection and preparation**

This assay may be used to measure Iodide.



It is the responsibility of the user to check the compatibility of the assay with the study matrix. To verify if there is no issue with the desired matrix, compare the data of the standard curve realised in water and in study matrix.

▷ **General precautions**

- ▶ All samples must be free from organic solvents prior to assay.
- ▶ Samples should be assayed immediately after collection or should be stored at -20°C .

▶ Reagent preparation

Each kit contains sufficient reagents for 96 wells. This allows for the construction of one standard curve in duplicate and the assay of 37 samples in duplicate.

All reagents need to be brought to room temperature (around +20°C) prior to the assay.

▶ Diluted nonradioactive iodide assay reagent A

The vial #D22076 has to be diluted four times with UltraPure water prior to use.

Use the following formula to determine the total volume for diluted nonradioactive iodide assay reagent A solution required:

$$(\text{Standards} + \text{Samples}) \times (\text{replicates}) \times 100\mu\text{L}$$

Then mix thoroughly by gentle inversion.

Stability: Use within the day

Example: If you want to assay the entire 96-well plate, you need 10mL of diluted nonradioactive iodide assay reagent A:

Dilute 2.5mL of #D22076 in 7.5mL of UltraPure water.

▷ Diluted nonradioactive iodide assay reagent B

Prepare the diluted nonradioactive iodide assay reagent B the same way as reagent A using the vial #D20076.

▷ Diluted nonradioactive iodide assay standards

Dilute 30 μ L of the vial #D06076 with 2970 μ L of UltraPure water; mix thoroughly by gentle inversion.

This 100x diluted solution is then diluted 10x a second time to create the stock solution: pipet 200 μ L of this solution with 1800 μ L of UltraPure water. Then mix thoroughly by gentle inversion.

The concentration of this stock solution is 2 μ M.

Prepare eight polypropylene tubes. Then prepare the standards by dilutions as follows:

| Standard | Volume of Standard stock solution at 2 μ M | Volume of UltraPure water | Standard concentration (nM) |
|----------|--|---------------------------|-----------------------------|
| S0 | - | 1mL | 0 |
| S1 | 350 μ L | 650 μ L | 700 |
| S2 | 300 μ L | 700 μ L | 600 |
| S3 | 250 μ L | 750 μ L | 500 |
| S4 | 200 μ L | 800 μ L | 400 |
| S5 | 150 μ L | 850 μ L | 300 |
| S6 | 100 μ L | 900 μ L | 200 |
| S7 | 50 μ L | 950 μ L | 100 |

Stability: Use within the day

▶ Assay procedure

It is recommended to perform the assays in duplicate following the instructions hereafter.

▶ Plate set-up

A plate set-up is suggested hereafter for the microplate procedure.

The contents of each well may be recorded on the template sheet provided at the end of this technical booklet.

| | 1 | 2 | 3 | 4 | 5 | 6 | 7 | 8 | 9 | 10 | 11 | 12 |
|---|----|----|----|---|---|---|---|---|---|----|----|----|
| A | Bk | S1 | S5 | * | * | * | * | * | * | * | * | * |
| B | Bk | S1 | S5 | * | * | * | * | * | * | * | * | * |
| C | Bk | S2 | S6 | * | * | * | * | * | * | * | * | * |
| D | Bk | S2 | S6 | * | * | * | * | * | * | * | * | * |
| E | S0 | S3 | S7 | * | * | * | * | * | * | * | * | * |
| F | S0 | S3 | S7 | * | * | * | * | * | * | * | * | * |
| G | S0 | S4 | * | * | * | * | * | * | * | * | * | * |
| H | S0 | S4 | * | * | * | * | * | * | * | * | * | * |

BK: Blank

S7-S0 : Standards

*: Samples

▷ Pipetting the reagents

All samples and reagents must reach room temperature prior to performing the assay.

Use different tips to pipet the buffers, nonradioactive iodide assay reagent A, nonradioactive iodide assay reagent B, standards and samples.

Before pipetting, equilibrate the pipette tips in each reagent. Do not touch the liquid already in the well when expelling with the pipette tip.

> **Blank**

Dispense 200 μL of water or matrix to prepare the standards to Blank (Bk) wells.

> **Nonradioactive iodide assay Standard**

Dispense 100 μL of each of the eight standards (S₀, then S₇ to S₁) in duplicate to appropriate wells. Start with the lowest concentration standard (S₀) and equilibrate the tip in the next higher standard before pipetting (S₇ to S₁).

> **Nonradioactive iodide assay Samples**

Dispense 100 μL in duplicate to appropriate wells.

> **Diluted nonradioactive iodide assay Reagent A**

Dispense 100 μL in each well except Blank (Bk) wells.

> **Diluted nonradioactive iodide assay Reagent B**

Dispense 100 μL in each well.

▷ Incubating the plate

Cover the plate with the cover sheet and incubate 30 minutes at +20°C in the dark.

▷ Reading the plate

- > Wipe the bottom of the plate with a paper towel, and make sure that no liquid has splashed outside the wells.
- > Read the plate at a wavelength between 414-420nm (yellow colour).

| Protocol in brief | | | | |
|---|-------|------------|----------|--------|
| Volume | Wells | Blank (BK) | Standard | Sample |
| Standard | | - | 100µL | - |
| Sample | | - | - | 100µL |
| Diluted reagent A | | - | 100µL | 100µL |
| Diluted reagent B | | 100µL | 100µL | 100µL |
| Water or Matrix used | | 200µL | - | - |
| Cover plate, incubate 30 min at 20 °C in dark | | | | |
| Read the plate at 414-420nm | | | | |

► Data analysis

Make sure that your plate reader has subtracted the absorbance readings of the blank well from the absorbance readings of the rest of the plate. If not, do it now.

- Calculate the average absorbance for each standard and sample.
- Calculate the log of each absorbance ($\log A$).
- For each standard, using a semi-log graph, plot the $\log A$ on y axis versus the concentration on x axis
- Use a linear regression.
- Calculate r^2 of your curve.
- To determine the concentration of your sample, find the absorbance value of each sample on the y axis.
- Read the corresponding value on the x axis which is the concentration of your unknown sample.
- Samples with a concentration greater than 700 nM should be re-assayed after dilution.

▶ **Acceptable range**

- > Regression coefficient $r^2 > 0.98$
- > Absorbance of S0 ≥ 0.800 A.U

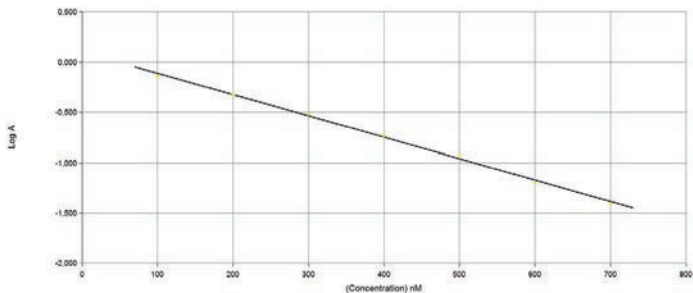
► Typical results

The following data are for demonstration purpose only. Your data may be different and still correct.

These data were obtained using all reagents as supplied in this kit under the following conditions: 30 minutes developing at +20°C, reading at 414 nm. A linear regression fitting was used to determine the concentrations..

| Standard | Expected concentration (nM) | Log A | Absorbance A.U |
|----------|-----------------------------|--------|----------------|
| S1 | 700 | -1.468 | 0.040 |
| S2 | 600 | -1.229 | 0.064 |
| S3 | 500 | -0.979 | 0.119 |
| S4 | 400 | -0.761 | 0.184 |
| S5 | 300 | -0.509 | 0.301 |
| S6 | 200 | -0.282 | 0.473 |
| S7 | 100 | -0.12 | 0.728 |

Typical Nonradioactive iodide standard curve



▶ Troubleshooting

- > **Absorbance values are too low:**
 - organic contamination of water,
 - one reagent has not been dispensed,
 - incorrect preparation / dilution,
 - assay performed before reagents reached room temperature,
 - incubation in wrong conditions (time or temperature).

- > **High signal and background in all wells:**
 - high ambient temperature.

- > **High dispersion of duplicates:**
 - poor pipetting technique.

These are a few examples of troubleshooting that may occur.

If you need further explanation, Bertin Pharma will be happy to assist you. Feel free to contact our technical support staff by phone (+33 (0)139 306 036), fax (+33 (0)139 306 299) or E-mail (bioreagent@bertinpharma.com), and be sure to indicate the batch number of the kit (see outside the box).

► Bibliography

Additional readings

List of publications quoting the use of this kit

1. Blazy P.
Iode
El-Aid JDID, 10 juin 2009
2. Role of Iodine in Metabolism.
*Chair and Department of Endocrinology, Jagiellonian University
Collegium Medicum 31-501 Krakow, 17 Kopernika St., Poland.*
3. Spitzweg C., Morris J. C.
The sodium iodide symporter: its pathophysiological and
therapeutic implications
22 october 2002
4. Porterfield S. P., Heindrich C.E.
The role of thyroid hormones in prenatal and neonatal
neurological development,
Endoc.Rev, 1993, pages 94-106
5. Anna Milanesi, Gregory A.Brent
Iodine and thyroid hormone synthesis, metabolism, and action.
*Molecular genetics and nutritional aspects of major and trace
minerals, 2017, pages 143-150*
6. Waltz F., Pillette L., Ambroise Y.
A nonradioactive iodide uptake assay for sodium iodide
symporter function.
Analytical Biochemistry, 16 Juillet 2009

7. Braun D., Schweizer I.
Thyroid Hormone transport and transporters,
Vitamins and hormones , 12 june 2017

| | | | | | | | | | | | | | | | | | |
|-----------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|
| 1 | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> |
| 2 | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> |
| 3 | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> |
| 4 | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> |
| 5 | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> |
| 6 | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> |
| 7 | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> |
| 8 | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> |
| 9 | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> |
| 10 | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> |
| 11 | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> |
| 12 | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> |
| A | | | | | | | | | | | | | | | | | |
| B | | | | | | | | | | | | | | | | | |
| C | | | | | | | | | | | | | | | | | |
| D | | | | | | | | | | | | | | | | | |
| E | | | | | | | | | | | | | | | | | |
| F | | | | | | | | | | | | | | | | | |
| G | | | | | | | | | | | | | | | | | |
| H | | | | | | | | | | | | | | | | | |



Bertin Pharma, over the last decades, has been developing and marketing over 100 biomarker assays, pre-analytical products, kits, antibodies and biochemicals thanks to its innovative work in research and development. Our core areas are orientated to inflammation, oxidative injury, endocrinology, diabetes, obesity, hypertension, neurodegenerative diseases, HIV, prion diseases, pharmacokinetics and metabolism.

Bertin Pharma is active worldwide either with direct sales or through our qualified and trained international distribution network from the United States to Japan.

We are able to provide you with local technical support to use at ease our products.

For further information, please send your request to
bioreagent@bertinpharma.com



care. innovate. deliver.

Parc d'activités du Pas du Lac - 10 bis avenue Ampère
F-78180 Montigny-le-Bretonneux - France
Tel: +33 (0)139 306 036 - Fax: +33 (0)139 306 299
bioreagent@bertinpharma.com - www.bertinpharma.com

—A **ENIM** Group Company—