

PRODUCT INFORMATION



Nitrite Standard

Item No. 780016

Purity: ≥98%

Supplied as: A lyophilized powder

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Reconstitute the lyophilized nitrite standard using 1.0 ml of assay buffer (20 mM potassium phosphate, pH 7.4). Be sure to mix in all powder that may cling to the vial stopper and vortex to ensure complete mixing. Dilute the standard 10-fold by adding 0.1 ml of the reconstituted standard to 0.9 ml of assay buffer. The concentration of the diluted solution is 200 μ M and is used for preparation of the nitrite standard curve as described in the table below. Store all stock solutions at 4°C; **do not freeze**. When stored under these conditions, the nitrite standard is stable for at least three months.

A typical standard curve can be prepared as follows:

Well (or Tube) Number	Nitrite Standard (μ l)	Assay Buffer (μ l)	Final Nitrite Concentration (μ M)
1	0	80	0
2	5	75	5
3	10	70	10
4	15	65	15
5	20	60	20
6	25	55	25
7	30	50	30
8	35	45	35

The concentration is calculated for a final volume of 200 μ l after the addition of 50 μ l each of Griess Reagent R1 and R2 (Item Nos. 780018 | 780020) as described in the kit booklet. When using a spectrophotometer for measuring absorbance values, the final volume and final concentrations may vary.

CAUTIONS

This product has been tested and formulated to work exclusively with the Nitrate/Nitrite Assay Kits from Cayman (Item Nos. 780001 | 780051). This product may not perform as described if used with other assay reagents or protocols.

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 03/07/2024

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM