

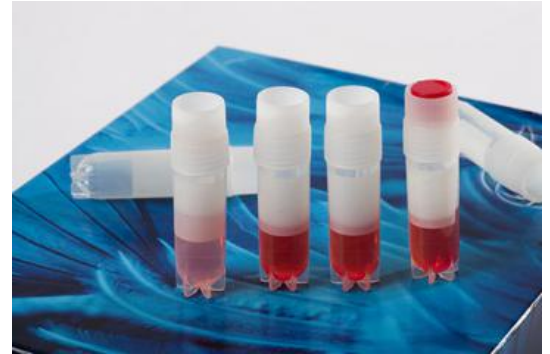


Sampling tubes with BHT and Indomethacin

Cat No: D31010 - 35 tubes

General Data

Shipping: room temp.



Application(s): BHT and Indomethacin sampling tubes are ready to use. Just collect your samples in classical blood collection tubes, and transfer 1 mL of sample per tube.

To be used to store samples after collection on a classical blood sampling tubes.

These tubes are not intended to be used to directly collect blood

Product Overview

These tubes are specially designed to be used for 1mL blood samples which will be assayed both for 8-isoprostanes and prostaglandins. See benefits of tubes D31007 and D31008.

The tubes are graduated, with a silicone washer sealed and external threads. They can be stored at temperatures as low as -80 degrees. Thus, BHT Indomethacin sampling tubes may be useful for any lipid peroxidation and COX metabolites studies.

FP/17/23

For research laboratory use only – Not for human diagnostic use.

Buyers agree to purchase the material subject to Purchasing Terms that can be found on our website. Seek appropriate training to safely handle this product under normal conditions. Use the recommended personal protective equipment to prevent chemical exposures.

Bertin Bioreagent does not make any other warranty or representation whatsoever whether expressed or implied, with respect to these products. In no event will Bertin Bioreagent be liable for incidental, consequential or punitive damages.

Contact Bertin Bioreagent

Parc d'Activités du Pas du Lac
10 bis avenue Ampère
78180 Montigny le Bretonneux - France
Tel.: +33 (0)139 306 036

<https://www.bertin-bioreagent.com/pa206/contact-us>
<https://www.bertin-bioreagent.com>