Product Information



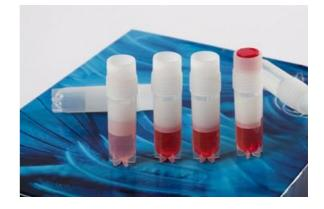


Sampling tubes with BHT

Cat No: D31007 - 35 tubes

General Data

Shipping:	room temp.
Stability:	Store at RT; shelf life 2 years maximum after production



Application(s): BHT 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

BHT Butylated hydroxytoluene (BHT) is a phenolic antioxidant and free radical scavenger. It has been shown to inhibit lipid peroxidation. Some biomarkers, like 8-isoprostanes, need to be collected on specific antioxidant in order to prevent artefactual formation or degradation of the biomarker. The BHT containing tubes have been designed to accept 1 mL of samples. 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 sampling tubes may be useful for any lipid peroxidation studies.

Scientific Literature

Antioxidant activity of vitamin E and related chain-breaking phenolic antioxidants in vitro", Journal of the American Chemical Society, 1981, volume 103

FP/27/17

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

Contact Bertin Pharma

10 bis avenue Ampère F-78180 Montigny-le-Bretonneux Tel.: +33 (0)139 306 036 bioreagent@bertinpharma.com **bioreagent.bertinpharma.com**