

Sampling tubes with PHMB

ORDERING INFORMATION

Catalogue Number: D31009

Size: 35 tubes or 100 tubes

Stability: 12 months

Storage: RT

Shipping: RT



- **CONCENTRATION:** 1 mM (final concentration)
- **SAMPLE SIZE:** 1 mL
- **APPLICATION:**

PHMB 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.
Those tubes are not intended to be used to directly collect blood.

- **DESCRIPTION:**

4-(hydroxymethyl)benzoic acid (PHMB) is used to inhibit cystein-protease and acts as a reversible thiol-specific blocker. While blocking the Acylated Ghrelin degradation, it allows the AcetylCholine Esterase AChE TM used in SPI-Bio EIA kits to work normally, in contrary to serine-protease inhibitors.

Neuropeptides and other short peptides are really sensitive to protease degradation. It is often necessary to add specific and potent inhibitor to avoid total loss of the biomarker. The PHMB containing tubes have been designed to accept 1 mL samples.

The cryotubes are graduated, with a silicone washer sealed and external threads. They can be stored at temperatures as low as -196°C but should be used only in the gas phase of liquid nitrogen.

Thus, PHMB sampling tubes may be useful to measure any biomarker cleaved by cystein-proteases.

- **ARTICLES:**

Boyer, P. D. (1954) J. Am. Chem. Soc. 76, 4331-4337,

For research laboratory use only.

Not for human diagnostic use.



Nov 09

Should you wish further information, do not hesitate to contact us.