

Product Information



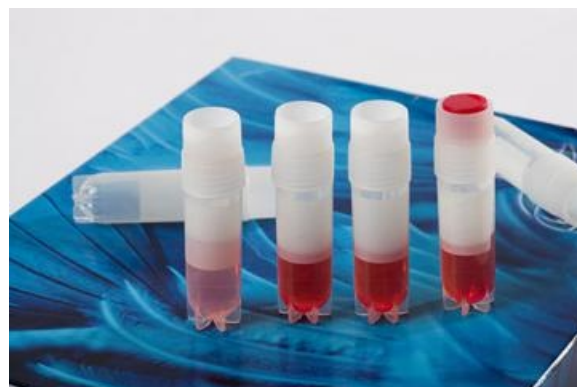
Sampling tubes with Indomethacin

Cat No: D31008 - 35 tubes

General Data

Shipping: room temp.

Stability: Store at RT; shelf life 2 years maximum after production



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

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

Product Overview

Indomethacin is a cyclooxygenase (COX) inhibitor that prevents ex-vivo Thromboxane B2 formation and other prostaglandins that can not be prevented using anticoagulant alone. The samples have to be transferred in Indomethacin sampling tubes as soon as possible after blood collection. For measuring some prostaglandins, it may be useful to block the ex-vivo formation of prostaglandin. The tubes are graduated, with a silicone washer sealed and external threads. They can be stored at temperatures as low as -80 degrees. Thus, indomethacin sampling tubes may be useful for any COX metabolites studies.

Scientific Literature

Ferreira S, Moncada S, Vane J (Jun 23 1971). "Indomethacin and aspirin abolish prostaglandin release from the spleen". Nat New Biol 231 (25): 237-9

FP/08/17

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