

Leukotriene B4

Leukotriene B4 is a pro-inflammatory dihydroxy fatty acid derived from arachidonic acid and produced from leukocytes in response to inflammation. Agonist of the BLT1 and BLT2 receptors. Promotes adhesion and activation of leukocytes at the endothelium, and stimulates production of pro-inflammatory cytokines and mediators. May be a factor in inflammatory diseases, and shows antiviral activity towards DNA viruses and retroviruses.

Materials Provided

Catalog Number:	92-1171
Quantity Supplied:	1 vial(s), 10 µg per vial

Description

DiscoverX control ligands have been functionally tested and validated for optimal use with all cell line targets and Assay Ready kits.

Product Information

Molecular Weight:	336.47 g/mol
Source:	Synthetic
Purity:	≥ 97%
Endotoxin Level:	N/A
Formulation:	N/A
Storage:	≤ -20°C Please avoid multiple freeze/thaw cycles.

Reagent Preparation

To avoid condensation, equilibrate the vial to ambient temperature before opening.

Stock Concentration:	300 µM
Reconstitution Volume:	0.100 mL
Reconstitution Solvent:	Ethanol

For Research Use Only

Ordering: +1.510.979.1415 option 4 or e-mail CustomerServiceDRX@eurofins.com
Technical support: +1.510.979.1415 option 5 or e-mail DRX_SupportUS@eurofinsUS.com
General product information: www.discoverx.com