

5-OxoETE

5-OxoETE is a keto acid analog of 5-hydroxyicosatetraenoic acid (5-HETE) that stimulates cytosolic calcium mobilization in neutrophils, actin polymerization, CD11b expression, and L-selectin shedding. It is a chemoattractant for eosinophils, neutrophils, basophils and monocytes, and an agonist of oxoeicosanoid (OXE) receptors (EC50=0.33nM).

Materials Provided

Catalog Number:	92-1042
Quantity Supplied:	1 vial(s), 25 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:	318.5 g/mol
Source:	Synthetic
Purity:	≥ 95%
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:	250 µM
Reconstitution Volume:	0.31390 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