

PAF

Platelet-activating factor (β -Acetyl- γ -O-alkyl-L- α -phosphatidylcholine, or PAF) is a phospholipid that mediates many leukocyte functions. Stimulates platelet aggregation and has roles in inflammation, anaphylaxis and wound repair. Produced by lymphocytes, platelets, and endothelial cells, and is also found to be secreted by tumor cells.

Materials Provided

Catalog Number:	92-1179
Quantity Supplied:	1 vial(s), 1 mg 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:	523.7 g/mol
Source:	Synthetic
Purity:	$\geq 99\%$
Endotoxin Level:	N/A
Formulation:	N/A
Storage:	$\leq -20^{\circ}\text{C}$ Please avoid multiple freeze/thaw cycles.

Reagent Preparation

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

Stock Concentration:	500 μM
Reconstitution Volume:	3.819 mL
Reconstitution Solvent:	DMSO

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