

Recombinant Human Complement C5a

Complement component 5a (C5a) is a pro-inflammatory protein fragment anaphylatoxin cleaved from complement component C5. It is a strong chemoattractant for leukocytes in stimulating an inflammatory response, and is recruited with the complement system network during an immune response to tissue damage or foreign objects. It is widely involved in many forms of inflammatory diseases, including asthma, ischemia-reperfusion injury, acute lung injury, and sepsis.

Materials Provided

Catalog Number:	92-1103
Quantity Supplied:	1 vial(s), 25 µg per vial
Reconstitution Buffer:	1 mL

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:	8300 g/mol
Source:	E. coli
Purity:	≥ 95% by SDS-PAGE
Endotoxin Level:	< 0.1 ng/µg (<1EU/µg).
Formulation:	Sterile filtered through a 0.2 micron filter. Lyophilized with no additives
Storage:	≤ -20°C Please avoid multiple freeze/thaw cycles.

Reagent Preparation

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

Stock Concentration:	25 µM
Reconstitution Volume:	0.120 mL
Reconstitution Solvent:	Reconstitution Buffer

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