

Anti-CD47 antibody

CD47 is an antiphagocytic ligand expressed by most cell types that binds the inhibitory receptor signal regulatory protein alpha (SIRP α) on phagocytic cells, sending a 'don't eat me signal' to the phagocyte.

Materials Provided

Catalog Number:	92-1341
Quantity Supplied:	1 vial(s), 20 μ g 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:	N/A
Source:	Mouse monoclonal antibody to human CD47; Isotype: IgG1/ κ
Purity:	N/A
Endotoxin Level:	N/A
Formulation:	200 μ g/ml of antibody purified from ascites fluid by Protein G chromatography. Prepared in
Storage:	2-8°C Please avoid multiple freeze/thaw cycles.

Reagent Preparation

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

Stock Concentration:	200 μ g/mL
Reconstitution Volume:	N/A
Reconstitution Solvent:	N/A

Handling Instructions:

stable for 24 months when stored at 2-8°C

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