

Apo-J Protein, Purified
Catalog Number: SIG-39085
Available Size: 50 uL

Description: Purified Apo-J / Clusterin Protein
(Formerly Signet Catalog Nos: 9085-005)

Intended Use: **Research Use Only (RUO)**

This product is sold for laboratory research use only, not for human or in-vivo use.

Form: Purified Protein (in PBS; no preservatives or carrier proteins)

Specificity: Purified mouse recombinant Apo-J/clusterin protein in PBS buffer. This product is provided at 1 mg/mL.

Suggested Working Dilution: The optimal working dilution should be determined for each specific assay condition.

Storage: Store at -20°C or below. Upon initial thawing, apportion into working aliquots and store at -20°C or below. Avoid repeated freeze-thaw cycles to prevent denaturing the antibody. Do not store in frost-free freezers.

Warranty/Conditions: Covance products may not be resold or modified for resale without prior written approval.

Product Revision Date: 4/4/2007