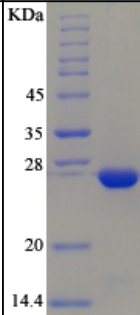


Recombinant Human APOBEC2 Protein Data Sheet

Catalog #	hRP-U0444-EF011
Size	25 µg
Protein Name	Homo sapiens apolipoprotein B mRNA editing enzyme, catalytic polypeptide-like 2
Protein Symbol	APOBEC2
Original Source	<i>Homo sapiens</i>
Expression System	<i>E.coli</i>
GenBank Accession #	NM_006789.1
Uniprot Accession #	Q9Y235
Description	Probable C to U editing enzyme whose physiological substrate is not yet known. Does not display detectable apoB mRNA editing. Has a low intrinsic cytidine deaminase activity. May play a role in the epigenetic regulation of gene expression through the process of active DNA demethylation.
Application	WB, ELISA, IP, antibody production, protein array
Fusion tag	N-His
Peptide Length	239aa(including fusion tag)
Molecular Weight	27.6kDa(including fusion tag)
pI	5.0
Activity	NA
Storage	Storage buffer: 20mM Tris.Cl, 50mM NaCl, 50% Glycerol, pH8.0. Store at -80°C and avoid repeated freeze-thaw cycles.
	
Reference:	• APOBEC2 is a monomer in solution: implications for APOBEC3G models • Excessive activity of apolipoprotein B mRNA editing enzyme catalytic polypeptide 2 (APOBEC2) contributes to liver and lung tumorigenesis. • An overview of cytidine deaminases

GeneCopoeia™
Expressway to Discovery

GeneCopoeia Inc.

9620 Medical Center Drive, Suite 101

Rockville, MD 20850

USA

Phone: 301-762-0888

Toll free: 1-301-762-0888

Fax: 301-762-3888

Web: www.genecopoeia.com

Inquiry: inquiry@genecopoeia.com

Technical Support: support@genecopoeia.com