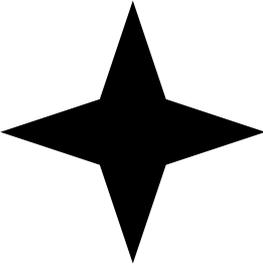


Recombinant Human ABHD14B Protein Data Sheet

Catalog #	hRP-V1020-EF011
Size	25 µg
Protein Name	Homo sapiens abhydrolase domain containing 14B
Protein Symbol	ABHD14B
Original Source	<i>Homo sapiens</i>
Expression System	<i>E.coli</i>
GenBank Accession #	NM_032750.1
Uniprot Accession #	Q96IU4
Description	Has hydrolase activity towards p-nitrophenyl butyrate (in vitro). May activate transcription.
Application	WB, ELISA, IP, antibody production, protein array
Fusion tag	N-His
Peptide Length	225aa
Molecular Weight	24.3kDa
pI	6.8
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:	Purification, crystallization and preliminary X-ray crystallographic analysis of human CCG1-interacting factor B.

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