Recombinant Human ASCL1 Protein Data Sheet

Catalog #	hRP-Z0740-EF012
Size	100 μg
Protein Name	Homo sapiens achaete-scute complex homolog 1 (Drosophila)
Protein Symbol	ASCL1
Original Source	Homo sapiens
Expression System	E.coli
GenBank Accession #	BC031299.1
Uniprot Accession #	P50553
Description	Transcriptional regulator. May play a role at early stages of development of specific neural lineages in most regions of the CNS, and of several lineages in the PNS. Essential for the generation of olfactory and autonomic neurons. Involved in the initiation of neuronal differentiation. Mediates transcription activation by binding to the E box (5'-CANNTG-3').
Application	WB, ELISA, IP, antibody production, protein array
Fusion tag	N-His
Peptide Length	251aa(including fusion tag)
Molecular Weight	27.4kDa(including fusion tag)
pI	9.0
Activity	NA
Storage	Storage buffer: 20mM Tris.Cl, 100mM NaCl, 50% Glycerol, pH8.0. Store at -80°C and avoid repeated freeze-thaw cycles.
	KDa 55 45 35 28



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

 $Technical \ Support: \underline{support@genecopoeia.com}$