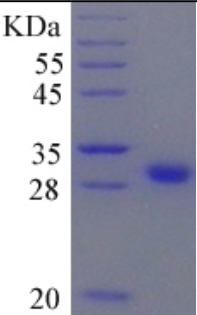


# Recombinant Human ASCL1 Protein Data Sheet

|                     |                                                                                                                                                                                                                                                                                                                                                                          |
|---------------------|--------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|
| Catalog #           | hRP-Z0740-EF012                                                                                                                                                                                                                                                                                                                                                          |
| Size                | 100 µg                                                                                                                                                                                                                                                                                                                                                                   |
| Protein Name        | Homo sapiens achaete-scute complex homolog 1 (Drosophila)                                                                                                                                                                                                                                                                                                                |
| Protein Symbol      | ASCL1                                                                                                                                                                                                                                                                                                                                                                    |
| Original Source     | <i>Homo sapiens</i>                                                                                                                                                                                                                                                                                                                                                      |
| Expression System   | <i>E.coli</i>                                                                                                                                                                                                                                                                                                                                                            |
| 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.                                                                                                                                                                                                                                                     |
|                     |                                                                                                                                                                                                                                                                                        |



## 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](http://www.genecopoeia.com)

Inquiry: [inquiry@genecopoeia.com](mailto:inquiry@genecopoeia.com)

Technical Support: [support@genecopoeia.com](mailto:support@genecopoeia.com)