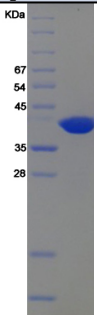


# Recombinant Human ADPRH Protein Data Sheet

|                     |                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                               |
|---------------------|---------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|
| Catalog #           | hRP-F0373-EF011                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                               |
| Size                | 25 µg                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                         |
| Protein Name        | [Protein ADP-ribosylarginine] hydrolase                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                       |
| Protein Symbol      | ADPRH                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                         |
| Original Source     | <i>Homo sapiens</i>                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                           |
| Expression System   | <i>E.coli</i>                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                 |
| GenBank Accession # | NM_001125.1                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                   |
| Uniprot Accession # | P54922                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                        |
| Description         | The enzyme encoded by this gene catalyzes removal of mono-ADP-ribose from arginine residues of proteins in the ADP-ribosylation cycle. Unlike the rat and mouse enzymes, which require DTT for maximal activity, the human enzyme is DTT-independent.                                                                                                                                                                                                                                                                                                                                                         |
| Application         | WB, ELISA, IP, antibody production, protein array                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                             |
| Fusion tag          | N-His                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                         |
| Peptide Length      | 372aa(including fusion tag)                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                   |
| Molecular Weight    | 41.4kDa(including fusion tag)                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                 |
| 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:          | <ul style="list-style-type: none"> <li>• <a href="#">Cloning and site-directed mutagenesis of human ADP-ribosylarginine hydrolase.</a></li> <li>• <a href="#">ADP-ribosylarginine hydrolases and ADP-ribosyltransferases. Partners in ADP-ribosylation cycles.</a></li> <li>• <a href="#">Crystal structure of dinitrogenase reductase-activating glycohydrolase (DraG) reveals conservation in the ADP-ribosylhydrolase fold and specific features in the ADP-ribose-binding pocket.</a></li> <li>• <a href="#">ADP-ribosylarginine hydrolase regulates cell proliferation and tumorigenesis.</a></li> </ul> |

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

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

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