

## SYVN1 (HRD1) Protein

Human recombinant protein expressed in Sf9 cells

**Catalog # S284-31G**

Lot # B1964-4

### Product Description

Recombinant human SYVN1 (HRD1) (226-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM\\_172230](#).

### Gene Aliases

SYVN1 (synoviolin); HRD1; DER3

### Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 25% glycerol.

### Storage and Stability

Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

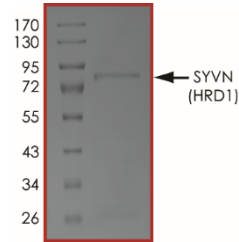
### Scientific Background

Synoviolin (SYVN1/HRD1) is a multispinning membrane protein with its C-terminal RING-H2 finger domain located in the cytoplasm. A member of the E3 ubiquitin-protein ligase family that accepts ubiquitin specifically from endoplasmic reticulum-associated UBC7 E2 ligase and transfers it to substrates, acting as a part of the ER-associated degradation (ERAD) system involved in ubiquitin-dependent degradation of misfolded endoplasmic reticulum proteins.

### References

1. Nadav E, et al: A novel mammalian endoplasmic reticulum ubiquitin ligase homologous to the yeast Hrd1. *Biochem Biophys Res Commun.* 2003 303(1):91-7.
2. Kikkert M, et al: Human HRD1 is an E3 ubiquitin ligase involved in degradation of proteins from the endoplasmic reticulum. *J Biol Chem.* 2004 279(5):3525-34.

### Purity



The purity of SYVN1 (HRD1) was determined to be **>80%** by densitometry. Approx. MW **~80kDa**.

## SYVN1 (HRD1) Protein

Human recombinant protein expressed in Sf9 cells

Catalog #	S284-31G
Lot #	B1964-4
Purity	>80%
Concentration	0.1 µg/µl
Stability	1yr at -70°C from date of shipment
Storage & Shipping	Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: [orders@signalchem.com](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

**FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.**