

## AMHR2 Protein

Recombinant human protein expressed in Sf9 cells

**Catalog # A38-35G**

Lot # A1243-3

### Product Description

Recombinant human AMHR2 (171-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. This gene accession number is [NM\\_020547](#).

### Gene Aliases

AMHR; MISR2; MISRII; MRII

### Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 50mM NaCl, 10mM glutathione, 0.1 mM 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.

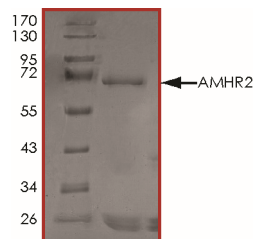
### Scientific Background

AMHR2 is a serine/threonine kinase with a single transmembrane domain belonging to the family of type II receptors for TGF-beta-related proteins which bind the ligand on their own but require the presence of a type I receptor for signal transduction. AMHR2 encodes the receptor for the anti-Mullerian hormone (AMH) which, in addition to testosterone which results in male sex differentiation and produced in the testes by different cells and has different effects (1). AMHR2 may be a regulator of neuronal pathways which vary in the male population and underlie subtle sex-linked biases (2).

### References

1. Imbeaud, S. et.al: Insensitivity to anti-mullerian hormone due to a mutation in the human anti-mullerian hormone receptor. *Nature Genet.* 11: 382-388, 1995.
2. Wang, P.-Y. et.al: Mullerian inhibiting substance contributes to sex-linked biases in the brain and behavior. *Proc. Nat. Acad. Sci.* 106: 7203-7208, 2009.

### Purity



The purity of AMHR2 protein was determined to be **>70%** by densitometry. Approx. MW **70 kDa**.

## AMHR2 Protein

Recombinant human protein expressed in Sf9 cells

Catalog #	A38-35G
Lot #	A1243-3
Purity	>70%
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.**