

IRAK3 Protein

Full length recombinant protein expressed in Sf9 cells

Catalog # 111-34G

Lot # V2250-6

Product Description

Recombinant full-length human IRAK3 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [BC057800](#).

Gene Aliases

IRAK-M

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 50mM 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.

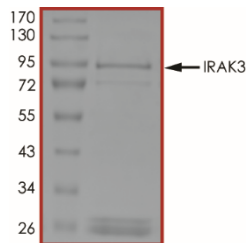
Scientific Background

IRAK3 or interleukin-1 receptor-associated kinase 3 is a member of the interleukin-1 receptor-associated kinase protein family which is essential components of the Toll/IL-R immune signal transduction pathways. IRAK3 is primarily expressed in monocytes and macrophages and functions as a negative regulator of Toll-like receptor signaling (1). IRAKM mediates critical aspects of innate immunity that result in an immunocompromised state during sepsis (2). Mutations in IRAK3 are associated with a susceptibility to asthma.

References

1. Del Fresno, C. et.al: Tumor cells deactivate human monocytes by up-regulating IL-1 receptor associated kinase-M expression via CD44 and TLR4. J. Immun. 174: 3032-3040, 2005.
2. Deng, J. C. et.al: Sepsis-induced suppression of lung innate immunity is mediated by IRAK-M. J. Clin. Invest. 116: 2532-2542, 2006.

Purity



The purity was determined to be **>70%** by densitometry. Approx. MW **96 kDa**.

IRAK3 Protein

Full length recombinant protein expressed in Sf9 cells

Catalog Number	111-34G
Specific Lot Number	V2250-6
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
www.signalchem.com

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