

HIF1B Protein

Full-length recombinant human protein expressed in Sf9 cells

Catalog # H07-30BG

Lot # Q2474-8

Product Description

Full-length recombinant human HIF1B was expressed by baculovirus in Sf9 cells using an N-terminal GST tag. The gene accession number is [NM_001668](#).

Gene Aliases

ARNT; bHLHe2; HIF-1-beta; HIF-1beta; HIF1-beta; HIF1BETA; TANGO

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.

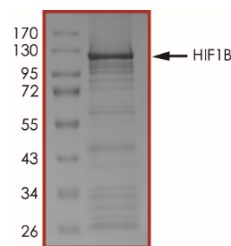
Scientific Background

The transcription factor Aryl hydrocarbon receptor nuclear translocator (ARNT), alternatively named Hypoxia-inducible factor (HIF)-1β, is part of the Hypoxia-inducible factor (HIF) pathway which mediates cellular adaptations to hypoxia and promotes tumour progression. Unlike HIF-1α that is regulated in an oxygen-dependent manner, HIF-1β is constitutively expressed. It has been shown that some tumour cells are capable to elevate HIF-1β in response to hypoxia to for a clonal advantage. HIF-1β binds HIF-1α, and forms the transcriptional active HIF-1 complex in the nucleus.

References

1. Mandl M, et al: (2015). The expression level of the transcription factor Aryl hydrocarbon receptor nuclear translocator (ARNT) determines cellular survival after radiation treatment. *Radiat Oncol*. 16;10:229.
2. Mandl M, et al (2013): Hypoxia-inducible factor-1β (HIF-1β) is upregulated in a HIF-1α-dependent manner in 518A2 human melanoma cells under hypoxic conditions. *Biochem Biophys Res Commun*. 26;434(1):166-72.

Purity



The purity was determined to be **>70%** by densitometry. Approx. MW **120kDa**.

HIF1B Protein

Full-length recombinant human protein expressed in Sf9 cells

Catalog #

H07-30BG

Lot #

Q2474-8

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.