

HIF-1 alpha antibody (mAb)

Catalog Nos: 61275, 61276

RRID: AB_2615045

Clone: ESEE122

Isotype: IgG1

Application(s): IHC, WB

Reactivity: Bovine, Human, Mouse, Rat

Quantities: 100 µg, 10 µg

Purification: Protein A Chromatography

Host: Mouse

Concentration: 1 µg/µl

Molecular Weight: 100 kDa

Background: HIF-1 (hypoxia inducible factor 1, alpha subunit, MOP1, PASD8, bHLHe78, PAS domain-containing protein 8) is a transcription factor found in mammalian cells cultured under reduced oxygen tension that plays an essential role in cellular and systemic homeostatic responses to hypoxia. HIF1 is a heterodimer composed of an alpha subunit and a beta subunit. The beta subunit has been identified as the aryl hydrocarbon receptor nuclear translocator (ARNT). This gene encodes the alpha subunit of HIF-1. Overexpression of a natural antisense transcript (aHIF) of this gene has been shown to be associated with nonpapillary renal carcinomas. Two alternative transcripts encoding different isoforms have been identified.

Immunogen: This HIF-1 alpha antibody was raised against a partial protein containing amino acid residues 329-530 of human HIF-1 alpha.

Buffer: Purified IgG in PBS with 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

This antibody is also available as an AbFlex[®] engineered recombinant antibody. For details on the corresponding AbFlex Recombinant Antibody, see Catalog No. 91139.

Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.