

TPX2 Protein

Full-length recombinant protein expressed in Sf9 cells

Catalog # T40-30H

Lot # N224-2

Product Description

Recombinant full-length human TPX2 protein was expressed by baculovirus in Sf9 insect cells using an N-terminal His tag. The gene accession number is [NM_012112](#). This product can be used for Aurora A activation in vitro.

Gene Aliases

DIL2, p100, DIL-2, HCTP4, FLS353, HCA519, REPP86, C20orf1, C20orf2, GD:C20orf1

Formulation

Recombinant protein stored in 50mM sodium phosphate, pH 7.0, 300mM NaCl, 150mM imidazole, 0.1mM PMSF, 0.25mM DTT, 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.

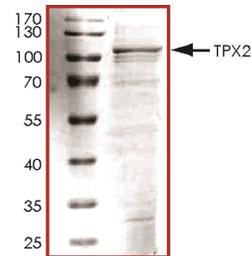
Scientific Background

TPX2 is a nuclear proliferation-associated protein whose expression is restricted to cell cycle phases S, G2, and M (1). It mainly expressed in the lung carcinoma cell lines and also found in adult placenta, skeletal muscle, thymus, testis, and small intestine and in fetal brain, liver, and kidney. TPX2 is required for targeting STK6 to the spindle apparatus, and that STK6 may regulate the function of TPX2 during spindle assembly. TPX2 is mainly function in mitotic spindle assembly likely contributes to its role in chromosome stability control and tumor suppression (2).

References

1. Heidebrecht, H. J. et.al: p100: a novel proliferation-associated nuclear protein specifically restricted to cell cycle phases S, G2, and M. Blood 90: 226-233, 1997.
2. Joukov, V.et.al: The BRCA1/BARD1 heterodimer modulates Ran-dependent mitotic spindle assembly. Cell 127: 539-552, 2006.

Purity



The purity of TPX2 was determined to be **>80%** by densitometry. Approx. MW **110kDa**.

TPX2 Protein

Full-length recombinant protein expressed in Sf9 cells

Catalog Number	T40-30H
Specific Lot Number	N224-2
Purity	>80%
Concentration	0.2 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.