

Importin β^{His}

D. melanogaster, Recombinant, *E. coli*

Cat. No.	Amount
PR-232	50 μg

Liquid. Supplied as a 1 mg/ml solution in 50 mM Tris-HCl, pH 7.5, 2 mM DTE, and 5 % glycerol.

Importins are involved in nuclear import and belong to a separate class of Ran-GTP binding proteins. Importin β binds the importin α -NLS complex or other adaptors to mediate the uptake of the substrate importin α/β complex through the nuclear pore complex. Importin α and β inhibit mitotic spindle assembly.

AVOID FREEZE/THAW CYCLES!

For in vitro use only!

weight: 98.69 kDa (884 amino acids)

Purity: > 95% by SDS-PAGE

Store: -80 °C

Selected References:

Dasso M. (2002) The Ran GTPase: Theme and Variations. *Current Biol.* **12**:502.

Macara I.G. (2001) Transport into and out the nucleus. *Microbiol. Biotech.* **65**:570.

Weis K. (1998) Importins and Exportins - How to get it in and out the nucleus. *Trends Biochem. Sci.* **23**:185.

Associated products available from Jena Bioscience

Non-hydrolyzable GTP-analogs (such as GTP γ S, GppNHp, NPE-caged-GTP...)
Molecular

Fluorescent GTPs (such as MANT, ANT, TNP, or GTP labeled with a dye of your choice)

Recombinant G-protein signaling reagents including GEFs, GAPs, small GTPases, heterotrimeric G-proteins, and many more...

For detailed information please view the sections on

<http://www.jenabioscience.com>