

Importin β^{His}

D. melanogaster, recombinant, *E. coli*

Cat. No.	Amount
PR-232	50 μg

For *in vitro* use only
Quality guaranteed for 12 months
Store at -80°C

Avoid freeze / thaw cycles

Form

Liquid. Supplied in 50 mM Tris-HCl pH 7.5, 2 mM DTE and 5 % glycerol.

Molecular Weight

98.7 kDa (884 amino acids)

Purity

> 95% by SDS-PAGE.

Description

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.

Selected References:

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

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

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

Associated products available from Jena Bioscience

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

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

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

For further information please visit www.jenabioscience.com.