

Importin α^{His} human, recombinant, *E. coli*

Cat. No.	Amount
PR-231	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, and 2 mM DTE.

Molecular Weight

57.86 kDa (529 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 α itself does not interact with the nuclear pore complex but functions as an adaptor that binds importin β . It is required to mediate binding of substrates to their receptors. Importin α binds directly to the nuclear localization signal of substrates.

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, small GTPases, heterotrimeric G-proteins, and many more...)

For further information please visit www.jenabioscience.com.