

## Importin $\alpha^{\text{His}}$ human, Recombinant, *E. coli*

Cat. No.	Amount
PR-231	50 $\mu\text{g}$

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

Importins are involved in nuclear import and belong to a separate class of Ran-GTP binding proteins. Importin  $\alpha$  itself does not interact with the nuclear pore complex but functions as an adaptor that binds importin  $\beta$ . It is required to mediate binding of substrates to their receptors. Importin  $\alpha$  binds directly to the nuclear localization signal of substrates.

AVOID FREEZE/THAW CYCLES!

### **For in vitro use only!**

**weight:** 57.86 kDa (529 amino acids)

**Purity:** > 95% by SDS-PAGE

**Store:** -80 °C

### **Selected References:**

Dasso M. (2002) The Ran GTPase: Theme and Variations. *Current Biology* **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 $\gamma$ S, GppCp, 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>