PICH (human), (recombinant)

PICH belongs to a family of SNF2-related nucleosome remodeling enzymes (DNA-dependent ATPases). These ATP-dependent motor proteins are major factors of chromatin structure and transcriptional control in vivo. SNF2-related factors have been implicated in a variety of cancers in humans.

Ordering Information

Order Online »

ENZ-PRT356-0050

50µl

Manuals, SDS & CofA

View Online »

Handling & Storage

Handling Thaw and freeze quickly, in room-temperature water bath and liquid nitrogen,

respectively; keep on ice when not in storage; split in small aliquots upon first thawing;

avoid excessive freeze-thaw cycles.

Long Term Storage -80°C

Shipping Dry Ice

Product Details

Alternative Name DNA excision repair protein ERCC-6-like, ERCC6L, PLK1-

ICH, RAD26L, Tumor antigen BJ-HCC-15

Formulation Liquid. In 20 mM Tris-HCl, pH 7.9, 150 mM NaCl, 15%

glycerol, 2 mM MgCl₂, 0.2 mM EDTA, 1 mM DTT, 10 mM beta-glycerophospate, 0.4 mM PMSF, 1 mM benzamidine

and 0.4 mg/ml recombinant human insulin.

MW ~143 kDa

Purity Detail Purified by immunoaffinity chromatography.

Source Produced in insect (Sf9) cells. Human PICH (aa 1-1250)

fused to 6xHis and FLAG-tag at the N-terminus.

UniProt ID Q2NKX8

Last modified: March 27, 2025



European Sales Office