## CSB (human), (recombinant)

CSB is an essential factor of trancription-coupled nucleotide excision repair (NER). It 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-PRT357-0050

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, Cockayne syndrome

B protein, ERCC6

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

glycerol, 2 mM MgCl<sub>2</sub>, 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** ~170 kDa

Purity Detail Purified by immunoaffinity chromatography.

**Source** Produced in insect (Sf9) cells. Human CSB (aa 1-1493)

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

UniProt ID Q03468-1

Last modified: March 27, 2025

