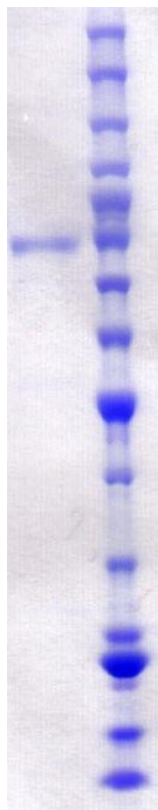


SUV420H2

Suppressor of Variegation 4-20 Homolog 2 (GST-His)

CATALOG NO.: HMT-21-349**LOT NO.:****DESCRIPTION:** Human recombinant SUV420H2 (residues 1-462; Genbank Accession # NM_032701.3; MW = 80.6 kDa) expressed as an N-terminal GST-His-fusion protein in *Sf9* insect cells.**PURITY:** >95% by SDS-PAGE**SUPPLIED AS:** $_ \mu\text{g}/\mu\text{L}$ in 50 mM Tris HCl, pH 7.5, 500 mM NaCl, 1 mM TCEP, 10% glycerol**STORAGE:** -70°C. Thaw quickly and store on ice before use. The remaining, unused, undiluted protein should be snap frozen, for example in a dry ice ethanol bath or liquid nitrogen. Minimize freeze/thaws if possible, but very low volume aliquots (<5 μl) or storage of diluted enzyme is not recommended.

Coomassie blue-stained SDS-PAGE (4-12% acrylamide) of 2 μg of RBC SUV420H2 (GST-His)
MW markers (right) are, from top, 220, 160, 120, 100, 90, 80, 70, 60, **50**, 40, 30, 25, **20**, 15, 10 kDa.

This product is not intended for therapeutic or diagnostic use in animals or in humans.

Reaction Biology1 Great Valley Parkway, Malvern PA, USA 19355
requests@reactionbiology.com www.reactionbiology.com