

S851-40N-30	30 µg
S851-40N-100	100 µg
S851-40N-1000	1 mg

Recombinant Human SHH

Recombinant human protein expressed in E. coli cells

Catalog # S851-40N

Lot # G3223-12

Product Description

Recombinant Human SHH was expressed in E. coli cells. The protein accession number is [Q15465](#)

Gene Aliases

SHH, HHG-1, Recombinant Human Sonic Hedgehog (SHH)

Endotoxin Level

<1.0 EU/µg of recombinant protein as determined by the LAL method

Formulation

Lyophilized from a 0.2 µm filtered solution in Phosphate buffer and NaCl

Reconstitution Protocol

A quick spin of the vial followed by reconstitution at 100 µg/mL in sterile PBS containing at least 0.1% human or bovine serum albumin.

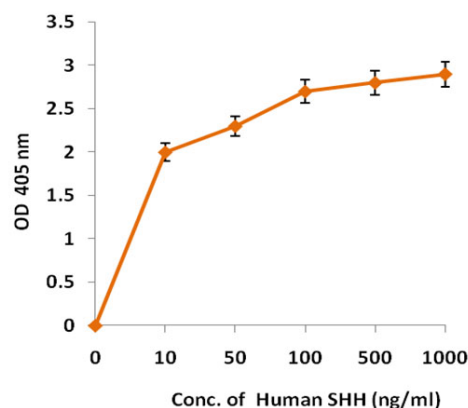
Storage and Stability

The lyophilized protein is stable for at least one year from date of receipt at -70°C. Upon reconstitution, this cytokine can be stored in working aliquots at 2° - 8°C for one month, or at -20°C for six months, with a carrier protein without detectable loss of activity. Avoid repeated freeze/thaw cycles.

Scientific Background

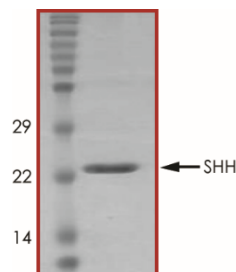
Highly conserved proteins, members of the Hedgehog family are widely represented throughout the animal kingdom. The three known mammalian Hedgehog proteins; Sonic (SHH), Desert (DHH) and Indian (IHH) hedgehog share a high degree of sequence and structural identity. SHH is specifically known to be expressed in fetal intestine, liver, lung and kidney. Recombinant Human SHH represents the N-product (N-terminal domain with Cys24 substituted by an Ile-Val-Ile sequence) which retains all the known signalling capabilities.

Specific Activity



Measured by its ability to induce alkaline phosphatase production by C3H10T1/2 mouse embryonic fibroblast cells. The ED50 for this effect is typically 0.1 - 0.4 µg/mL.

Purity



Recombinant Human SHH resolved on a 17% SDS-PAGE gel under reducing conditions and stained with Coomassie Brilliant Blue G-250

Approx. MW **20.1 kDa**

Recombinant Human SHH

Recombinant human protein expressed in E. coli cells

Catalog Number	S851-40N
Specific Lot Number	G3223-12
Purity	95%
Stability	6 mos. at -20°C from date of shipment
Storage & Shipping	The lyophilized protein is stable for at least one year from date of receipt at -70°C. Upon reconstitution, this cytokine can be stored in working aliquots at 2° - 8°C for one month, or at -20°C for six months, with a carrier protein without detectable loss of activity. Avoid repeated freeze/thaw cycles.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.