

Catalog #	Aliquot Size
D834-40N-50	50 μ a
D834-40N-250	250 μ g
D834-40N-1000	1 mg

Recombinant Human DHH

Recombinant human protein expressed in E. coli cells

Catalog # D834-40N

Lot # G3211-25

Product Description

Recombinant Human DHH was expressed in E. coli cells. The protein accession number is [O43323](#)

Gene Aliases

Desert Hedgehog Protein, DHH
Recombinant Human Desert Hedgehog (DHH)

Endotoxin Level

<1.0 EU/ μ g of recombinant protein as determined by the LAL method.

Formulation

Lyophilized from a 0.2 μ m filtered solution in PBS with 0.2 mM DTT

Reconstitution Protocol

Reconstitute at 500 μ g/mL in sterile PBS

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.

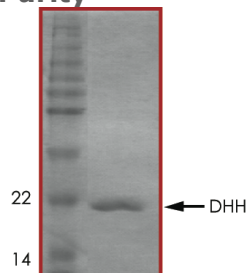
Scientific Background

The Desert Hedgehog protein is one of the three Hedgehog family proteins that is highly conserved and plays a role in many developmental processes in animals. DHH become biologically active following its autocatalytic cleavage. There is a high degree of residue homology between Sonic Hedgehog (SHH), Indian Hedgehog (IHH), and DHH. Despite their individual levels of expression and unique biological roles, the Hedgehog family members all use the same signalling pathway, and therefore, may be used interchangeably for some experimental purposes.

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 15-45 μ g/ml

Purity



Recombinant Human DHH resolved on a 15% SDS-PAGE gel under reducing conditions and stained with Coomassie Brilliant Blue G-250

Approx. MW **20.0 kDa**

Recombinant Human DHH

Recombinant human protein expressed in E. coli cells

Catalog Number	D834-40N
Specific Lot Number	G3211-25
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.