

Echelon Biosciences Inc.

Biotinylated Versican G1 Domain (bVG1)

G-HA02

Support: echelon@echelon-inc.com

Description:

Biotinylated N-terminal versican G1 domain.

Properties:

Size – 50 µg

Form – Liquid

Storage Buffer – PBS (0.01 M phosphate buffer, 0.0027 M potassium chloride and 0.137 M sodium chloride, pH 7.4).

Source – Recombinant human protein expressed in E.coli¹

Concentration – Please see Certificate of Analysis for lot specific information

Purity – Please see Certificate of Analysis for lot specific information

Specificity – No significant cross reactivity to other glycosaminoglycans based on data from a direct ELISA

Storage – Store at -80 °C. Protein is stable for at least 6 months from date of shipment. Avoid freeze thaw cycles.

Applications:

ICC – 5 µg/mL

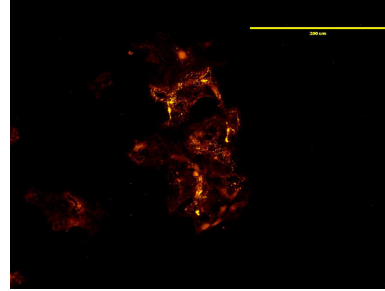
ELISA – 0.05-1.25 µg/mL

Other in vitro and cellular applications are possible using this protein but have not been verified by Echelon Biosciences. Optimal usage concentration should be optimized by researcher.

Background:

Versican is one of the many hyaluronan binding proteins that belong to the link module superfamily. Versican binds hyaluronan through the specific binding domain called “link module”. The G1 domain of versican contains an immunoglobulin domain and two contiguous link modules.

Data: ICC



HEK293 cells stained with bVG1 and fluorophore-conjugated Streptavidin.
Scale bar = 200 µm

References:

1. Seyfried, N.T., et al., Expression and purification of functionally active hyaluronan-binding domains from human cartilage link protein, aggrecan and versican: formation of ternary complexes with defined hyaluronan oligosaccharides. *J Biol Chem*, 2005. 280(7): p. 5435-48.
2. Rivas, F., et al., Label-free analysis of physiological hyaluronan size distribution with a solid-state nanopore sensor. *Nat Commun*, 2018. 9(1): p. 1037.
3. Yuan, H., et al., Determination of hyaluronan molecular mass distribution in human breast milk. *Anal Biochem*, 2015. 474: p. 78-88.

Related Products:

Product	Catalog Number
Assays	
HA Quantification ELISAs	K-1200, K-4800
Hyaluronidase Activity ELISA	K-6000
Compounds	
BODIPY-HA	H-025F, H-250F, H-700F
Texas Red-HA	H-025R, H-250R, H-700R
Select-HA™	HYA-50kEF-1, HYA-500kEF-1, HYA-601kEF-1, HYA-1000kEF-1
Biotin Select-HA™	HYA-B50-200, HYA-B250-200, HYA-B500-200, HYA-B100-200
Select-HA Ladder™	HYA-MGLAD-20, HYA-HILAD-20, HYA-LOLAD-20
nanoHA™	HYA-NAN05-1
HAase Inhibitor	B-0601
HA Synthese Inhibitor	JK201388
HA Binding Proteins	
Versican G1 Domain	G-HA01, G-HA03

Technical Data Sheet Rev. 3, 05-20-20 - For research use only. Not intended or approved for diagnostic or therapeutic use.