

Histidine Tagged Versican G1 Domain (His-VG1)

G-HA03

Support: echelon@echelon-inc.com

Description:

C-terminal 12X Histidine tagged Versican G1 domain.

Applications:

ELISA – 25 to 200 ng/mL

Other in vitro and cellular applications are possible using this antibody but have not been verified by Echelon Biosciences.

Properties:

Size – 50 µg

Form – Liquid

Storage Buffer – PBS (10 mM phosphate buffer, 2.7 mM potassium chloride and 137 mM sodium chloride, pH 7.4).

Source – 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.

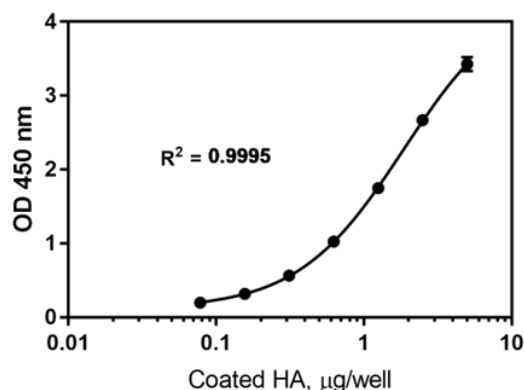
Specificity:

The Versican G1 Domain binds to hyaluronan (HA). This makes it a valuable tool for detecting hyaluronan in various biological samples.

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: Direct ELISA



HA coated on 96-well plate detected by 25 ng/mL his-VG1.

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.

Related Products:

Products	Catalog Number
Assays	
HA Quantification Assays	K-1200, K-4800, K-5800
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
HAse / HA Synthase Inhibitor	B-0601 / JK201388
HA Binding Proteins	
Versican G1 Domain	G-HA01, G-HA02