



Echelon Biosciences Inc.  
675 Arapeen Drive, Suite 302  
Salt Lake City, UT 84108  
Telephone 866-588-0455  
Fax 801-588-0497  
echelon@echelon-inc.com  
www.echelon-inc.com

## Technical Data Sheet

**For research use only**  
Not intended or approved for  
diagnostic or therapeutic use.

**Product Name:** Versican G1 Domain (VG1)

**Product Number:** G-HA01

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

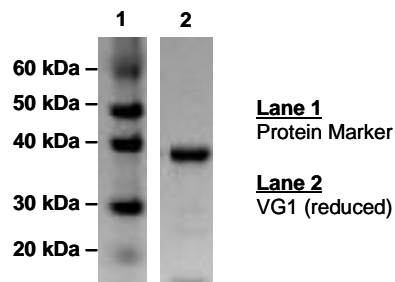
**Description:** Versican is one of the many hyaluronan binding proteins that belong to the link module superfamily. Versican binds hyaluronan through the specific binding domain called “link module”, in which, it contains an immunoglobulin domain and two contiguous link modules. Echelon VG1 contains only the G1 domain from versican.

**Source:** Recombinant human versican G1 domain expressed in *E.coli*.

**Concentration:** 50 µL at 1 mg/mL in PBS (0.01 M phosphate buffer, 0.0027 M potassium chloride and 0.137 M sodium chloride, pH 7.4).

**Suggested Use:** 0.5-1 µg/mL in ELISA. Optimal usage concentration should be determined by researcher. *Other applications have not been tested.*

**Purity:** On SDS-PAGE commassie staining gel under reducing condition, the VG1 showed a band at ~36.7 kDa with purity ≥ 80%.



**References:** Seyfried NT, McVey GF, Almond A, Mahoney DJ, Dudhia J, Day AJ. 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 Feb 18;280(7):5435-48.

**Related Products:** Hyaluronan Enzyme-Linked Immunosorbent Assay (K-1200)  
Hyaluronic Acid (HA) Screening Service (T-1200)  
Collagen IV ELISA Kit (K-1700)  
Hyaluronic Acid Sandwich ELISA (K-4800)