

GroPep Bioreagents

Bovine Betacellulin antiserum (Rabbit)

Description

Bovine betacellulin antiserum was prepared by immunising rabbits with recombinant bovine betacellulin. It has 100% cross-reactivity with bovine betacellulin but not to betacellulin from other species. It does not recognise human epidermal growth factor (EGF) or human transforming growth factor- α (TGF- α).

References:	Bastian SEP <i>et al</i> (2001) Journal of Endocrinology 163 , 203-212
Immunogen:	Bovine betacellulin
Source:	Rabbit antiserum
Specificity:	Cross-reacts 100% with bovine betacellulin, <1% with rat betacellulin. There was no cross-reactivity to human EGF or TGF- α .
Appearance:	Lyophilized powder
Storage/Stability:	At least 2 years at 2 - 4°C (lyophilized). After reconstitution store at -20°C or -80°C Avoid freeze-thaw cycles
Reconstitution:	Dissolve the vial contents in 500 μ l sterile water and stand for 10 minutes at room temperature. Dilute further in a buffer suitable for the application.
Application and Titre:	Each application and titre should be determined in house but as a guideline: Radioimmunoassay: 1:300,000 Western blotting: 1:1000

Product Codes and Pricing

Bovine betacellulin antiserum **100 μ l** **PAAR1**

Related Products: Bovine betacellulin
Rat betacellulin antiserum (Rabbit)

****NOT FOR USE IN HUMANS****

GroPep Bioreagents Pty Ltd
51 West Thebarton Road
Thebarton SA 5031
Australia

ABN 93 147 032 166

Telephone: +61 8 7222 1051

Postal Address:
PO Box 10065
Adelaide Business Centre
SA 5000
Australia

Email:
info@gropep.com

Internet:
www.gropep.com