



GroPep Bioreagents



Rat Betacellulin antiserum (Rabbit)

Description

Rat betacellulin antiserum was prepared by immunising rabbits with a 16 amino acid peptide derived from a sequence downstream from the N terminus of mature rat betacellulin. It has 100% cross-reactivity with rat betacellulin but not to bovine or human betacellulin. It does not recognise human epidermal growth factor (EGF) or human transforming growth factor- α (TGF- α).

References:

Dunbar AJ & Goddard C (2000) International Journal of Biochemistry and Cell Biology **32**, 805-815

Immunogen:

Derived from UniProtKB/Swiss-Prot sequence ID Q9JIM4 rat betacellulin. The peptide sequence is amino-acids 50-65 (rat betacellulin₁₉₋₃₄) conjugated to diphtheria toxoid.

Source:

Rabbit antiserum

Specificity:

Cross-reacts 100% with rat betacellulin, <1% with bovine 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 100 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:

Western blotting: 1:1000

Product Codes and Pricing

Rat betacellulin antiserum 100 μ l PAAQ1

Related Products:

Bovine betacellulin
Bovine 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