

PRODUCT INFORMATION



Guanylate Cyclase β_1 subunit (soluble) Blocking Peptide

Item No. 360897

Overview and Properties

Contents:	A <i>batch specific</i> quantity of peptide, consisting of the rat soluble guanylate cyclase β_1 subunit. Batch specific information can be found online or by contacting Technical Support.
Synonyms:	sGC β_1 -subunit
Immunogen:	Synthetic peptide from the N-terminal 250 amino acids {rat}
Uniprot No.:	P20595
Form:	Solid
Storage:	-20°C (as supplied)
Storage Buffer:	Re-suspend peptide in 200 μ l ddH ₂ O
Stability:	\geq 3 years (as supplied)
Purity:	\geq 95%

Applications

This blocking peptide can be used in conjunction with Cayman's Guanylate cyclase β_1 subunit (soluble) Polyclonal Antibody (Item No. 160897) to block protein-antibody complex formation during immunochemical detection of guanylate cyclase β_1 subunit.

Procedure:

1. Mix 50 μ l Guanylate Cyclase β_1 subunit Polyclonal Antibody (Item No. 160897) and 20 μ l blocking peptide together in a microfuge tube.*
2. Incubate for one hour at room temperature with occasional mixing prior to further dilution and application of the mixture to the immunoblot.
3. Dilute the mixture to the final working antibody concentration and apply to the slide or membrane as usual.

*This is a recommended mixture. The minimum amount of peptide needed for complete blocking has not been precisely determined and may vary depending on the sample being analyzed. The amount of peptide required may need to be increased if sufficient blocking does not occur

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/15/2023

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM