

PRODUCT INFORMATION



5-OxoETE Receptor Blocking Peptide

Item No. 10006618

Overview and Properties

| | |
|------------------------|--|
| Contents: | This vial contains 200 µg of peptide. |
| Synonyms: | R527, TG1019 |
| Form: | Liquid |
| Storage: | -20°C (as supplied) |
| Storage Buffer: | 200 µl TBS, pH 7.4, with 0.1% BSA and 0.02% sodium azide |
| Stability: | ≥3 years |

Procedures

This vial contains 200 µg peptide in 200 µl TBS, pH 7.4, with 0.1% BSA and 0.02% sodium azide. The peptide is derived from the human 5-OxoETE receptor C-terminal amino acids 408-423 (KVQGEVSLEKEGSSQG). This blocking peptide can be used in conjunction with Cayman's 5-OxoETE Receptor Polyclonal Antibody (Item No. 100025) to block protein-antibody complex formation during immunochemical analysis of 5-OxoETE receptor.

To block antibody/protein complex formation, the following procedure is recommended:

1. Mix the 5-OxoETE Receptor Polyclonal Antibody (Item No. 100025) and blocking peptide together in a 2:1 (v/v) ratio in a microfuge tube. For example, mix 20 µl of antibody and 20 µl of peptide.*
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, 10/16/2023

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