

PRODUCT INFORMATION



CysLT₂ Receptor (C-Term) Blocking Peptide Item No. 320550

Overview and Properties

Contents:	This vial contains 200 µg of peptide.
Synonym:	Cysteinyl Leukotriene Receptor 2
Storage:	-20°C (as supplied)
Storage Buffer:	200 µl TBS, pH 7.4, containing 0.1% BSA and 0.02% sodium azide
Stability:	≥3 years

Description

The cysteinyl leukotrienes (cysLTs; LTC₄, LTD₄, and LTE₄) contract airway and pulmonary vascular smooth muscle, increase vascular permeability, and stimulate mucus secretion, thereby playing a major role in asthma.¹⁻⁴ LTC₄, LTD₄, and LTE₄ mediate their actions *via* at least two receptors designated CysLT₁ and CysLT₂.¹ Cloning of the human CysLT₂ receptor reveals it is a 346 amino acid protein with 38% homology to the CysLT₁ receptor.^{5,6} The rank order of binding for LTs to the cloned receptor, as determined using a radioligand binding assay, is LTC₄ = LTD₄ >> LTE₄.⁵ The mRNA for the human CysLT₂ receptor is expressed in lung macrophages, airway smooth muscle, cardiac Purkinje cells, adrenal medulla cells, peripheral blood leukocytes, spleen, placenta, and brain.^{5,7}

Procedures

This vial contains 200 µg peptide in 200 µl TBS, pH 7.4, containing 0.1% BSA and 0.02% sodium azide. The CysLT₂ receptor (C-term) blocking peptide (human CysLT₂ receptor amino acids 330-346) can be used in conjunction with Cayman's CysLT₂ Receptor (C-Term) Polyclonal Antibody (Item No. 120550) to block protein-antibody complex formation during immunochemical analysis of the CysLT₂ receptor.

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

1. Mix the CysLT₂ Receptor (C-Term) Polyclonal Antibody (Item No. 120550) and blocking peptide together in a 1:1 (v/v) ratio in a microfuge tube. For example, mix 40 µl of antibody and 40 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.

References

1. Gorenne, I., Norel, X., and Brink, C. *Trends Pharmacol. Sci.* **17**, 342-343 (1996).
2. Dahlén, S.-E., Hansson, G., Hedqvist, P., et al. *Proc. Natl. Acad. Sci. USA* **80**, 1712-1716 (1983).
3. Busse, W.W. *Clin. Exp. Allergy* **26**, 868-879 (1996).
4. Hedqvist, P., Dahlén, S.-E., Gustafsson, L., et al. *Acta Physiol. Scand.* **110**, 331-333 (1980).
5. Heise, C.E., O'Dowd, B.F., Figueroa, D.J., et al. *J. Biol. Chem.* **275**, 30531-30536 (2000).
6. Lynch, K.R., O'Neill, G.P., Liu, Q., et al. *Nature* **399**, 789-793 (1999).
7. Takasaki, J., Kamohara, M., Matsumoto, M., et al. *Biochem. Biophys. Res. Commun.* **274**, 316-322 (2000).

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, 12/14/2023

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