

CERTIFICATE of ANALYSIS



BLT1 Receptor Monoclonal Antibody (Clone 7B1)

Item No. 120111 • Batch No. 0630233 • Expires 24JAN2028

Overview

Formulation:	200 µg of Protein-G purified IgG
Storage/stability:	-20°C/1yr
Applications:	FC, IF, and IHC of frozen sections
Cross reactivity:	(+)BLT1 receptor(-)BLT2 receptor(-)CysLT1 receptor(-)CysLT2 receptor
Species reactivity:	(+)Human BLT1 receptor

Reviewed and approved by: Kyle Bale

Approved date: 03NOV2021

A handwritten signature in black ink that reads "Kyle Bale".

WARNING

THIS PRODUCT IS FOR RESEARCH USE - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE. IT IS THE RESPONSIBILITY OF THE PURCHASER TO DETERMINE SUITABILITY FOR OTHER APPLICATIONS.

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, 10/12/2018

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