

# CERTIFICATE of ANALYSIS



## BLT1 Receptor Polyclonal Antiserum

Item No. 100019 • Batch No. 0431342 • Expires 17JUN2027

### Overview

**Formulation:** 100 µl lyophilized antiserum  
**Storage/Stability:** -20°C/1yr  
**Applications:** CF, FC, IHC (formalin-fixed paraffin-embedded tissue), and WB  
**Species reactivity:** (+)Bovine, (+)Human, (-)Mouse

Reviewed and approved by: Paul Domanski

Approved date: 07SEP2016

A handwritten signature in black ink that reads "Paul Domanski".

**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