

CERTIFICATE of ANALYSIS



5-Lipoxygenase (Phospho-Ser523) Polyclonal Antibody

Item No. 10007820 • Batch No. 0679053 • Expires 17MAY2024

Overview

Formulation:	Affinity-purified rabbit polyclonal antibody; stored in 10 mM HEPES (pH 7.5), 150 mM NaCl, 100 µg per ml BSA and 50% glycerol
Storage/Stability:	-20°C/1yr
Applications:	WB
Species reactivity:	(+)Human 5-<wbr/>LO, (+)Non-<wbr/>human primate 5-<wbr/>LO, (+)Rat 5-<wbr/>LO

Reviewed and approved by: Judy Hines

Approved date: 17MAY2023

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