

CERTIFICATE of ANALYSIS



Citrullinated Collagen Type II Chimeric Rabbit Monoclonal Antibody (Clone ACC1)

Item No. 39297 • Batch No. 0702625 • Expires 15FEB2025

Overview

Formulation:	200 µg of protein A-affinity purified monoclonal antibody PBS with 0.02% Proclin 300.
Concentration:	1 mg/ml
Storage/Stability:	-20°C/1yr
Applications:	ELISA, IHC
Cross reactivity:	(+)Citrullinated form of the triple helical collagen type II C1 epitope (amino acids 359-369), (+)Denatured collagen type II, (+)Denatured collagen type II triple helical C1 epitope (amino acids 359-369)
Species reactivity:	(+)Bovine, (+)Human, (+)Mouse

Reviewed and approved by: Emily Tucker

Approved date: 01MAR2024

A handwritten signature in black ink, appearing to read "Emily Tucker".

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