

PRODUCT INFORMATION



Normal Goat Serum

Item No. 10006577

Overview and Properties

Contents:	This vial contains 20 ml of 10% normal goat serum in PBS, pH 7.4, with 0.04% sodium azide.
Storage:	-20°C
Stability:	≥3 years
Applications:	This solution can be used (without further dilution) as a blocking agent for immunocytochemistry or immunohistochemistry when using a secondary antibody generated in goat. For example, incubating a sample with this normal serum would be applicable prior to use of Cayman's Goat Anti-Rabbit IgG FITC (Item No. 10006588).

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, 10/16/2023

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