

Product Information



Transcription Factor Goat Anti-Mouse HRP Conjugate

Catalog No. 10009279

Contents: This vial contains 100 µl of HRP-conjugated Goat Anti-Mouse IgG in 50% glycerol.

Stability: ≥1 year at -20°C

Laboratory Procedures

This antibody has been specifically formulated to be used in Transcription Factor Assays. Prepare the HRP-conjugated Goat Anti-Mouse HRP Conjugate by diluting 1:100 in 1X Antibody Binding Buffer (ABB) as specified in the table below. Add 100 µl of diluted Goat Anti-Mouse HRP Conjugate at the appropriate stage in the assay to each well of the supplied 96-well plate except the Blank (Blk) wells. Discard any unused antibody.

Component	Volume/Well	Volume/Strip	Volume/96-Well Plate
UltraPure water	89 µl	712 µl	8,544 µl
10X Antibody Binding Buffer	10 µl	80 µl	960 µl
Transcription Factor p53 Primary Antibody	1 µl	8 µl	96 µl
Total required	100 µl	800 µl	9,600 µl

CAUTIONS

For laboratory research use only. This reagent is not for human or veterinary use. This product has been tested and formulated to work exclusively with kits from Cayman.

Related Products

96-Well Cover Sheets - Cat. No. 400012 • Tween 20 - Cat. No. 400035 • Wash Buffer Concentrate (400X) - Cat. No. 400062 • p53 Transcription Factor Assay Kit - Cat. No. 600020 • Transcription Factor p53 96-Well Plate - Cat. No. 600021 • Transcription Factor p53 Positive Control - Cat. No. 600022 • Transcription Factor p53 Primary Antibody - Cat. No. 600023 • Transcription Factor p53 Specific Competitor dsDNA - Cat. No. 600024 • Transcription Factor Binding Assay Buffer (4X) - Cat. No. 10006880 • Transcription Factor Antibody Binding Buffer (10X) - Cat. No. 10006882 • Transcription Factor Developing Solution - Cat. No. 10006888 • Transcription Factor Stop Solution - Cat. No. 10006889 • Transcription Factor Reagent A - Cat. No. 10007472

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY; NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent under separate cover to the MSDS supervisor at your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees. Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy located on our website and in our catalog.**

Copyright Cayman Chemical Company, 10/21/2008