

Product Information



Transcription Factor p53 Primary Antibody

Item No. 600023

Synonym: TP53, TRP53, LFS1
Host: Mouse
Applications: p53 Transcription Factor Assay Kits (Item Nos. 600020 and 600030)
Stability: ≥1 year at -20°C

Laboratory Procedures

This antibody has been specifically formulated to be used in Cayman's p53 Transcription Factor Assay Kits (Item Nos. 600020 and 600030). Prepare the primary antibody by diluting 1:100 in 1X Transcription Factor Antibody Binding Buffer diluted according to the procedure as outlined below. Add 100 µl of diluted p53 Antibody to each well except the Blank (Blk) wells. Discard any unused diluted primary 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 the p53 Transcription Factor Assay Kits from Cayman (Item Nos. 600020 and 600030). This product may not perform as described if used with other assay reagents or protocols.

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/600023

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.

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 Safety Data Sheet, which has been sent via email to 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, 06/14/2013