

Product Information



Transcription Factor p53 Specific Competitor dsDNA

Item No. 600024

- Contents:** This vial contains 120 µl specific competitor double stranded oligonucleotide
- Stability:** ≥1 year at -20°C.
- Storage:** It is recommended that the p53 specific competitor dsDNA be stored at -20°C.
- Application:** The provided competitor double stranded oligonucleotide is a specific competitor for p53 transcription factor binding and demonstrates specificity of the assay. This competitor dsDNA competes with the sequence bound to the plate for p53 binding. The competitor oligonucleotide is not biotinylated and therefore will not bind to the plate. In the control nuclear extracts where p53 specific competitor oligonucleotide has been added, the signal intensity is significantly reduced. The competitor does not need to be run with every assay.

Laboratory Procedures

Add 10 µl of the p53 specific competitor dsDNA to each well as specified in the p53 Transcription Factor Assay Kit booklets (Item Nos. 600020 and 600030). These wells should show reduced signal intensity when compared to positive control wells where competitor dsDNA has not been added.

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 Assays 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/600024

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