

Product Information



TRF2 Polyclonal Antibody

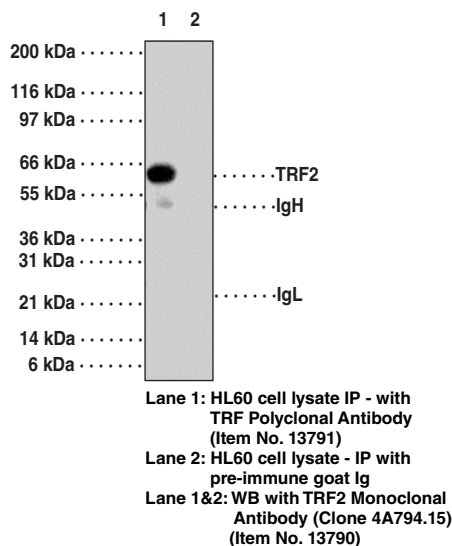
Item No. 13791

| | |
|--------------------------|--|
| Contents: | This vial contains 100 µg of Protein G purified IgG in 200 µl PBS containing 0.05% BSA and 0.05% sodium azide. |
| Synonyms: | Telomeric Repeat Binding Factor 2, TERF2, TRBF2, TTAGGG Repeat Binding Factor 2 |
| Antigen: | Full length human TRF2 |
| GenBank | |
| Accession No.: | NP_005643 |
| Host: | Goat |
| Cross Reactivity: | (+) Human and mouse TRF2 |
| Storage: | ≤6 months at 4°C; ≥1 year at -20°C |
| Applications: | ChIP, immunocytochemistry, immunofluorescence, immunohistochemistry (IHC), immunoprecipitation (IP), and western blot (WB); the recommended starting concentration for IHC is 15 µg/ml, 2 µg/1x10 ⁶ cells for IP, and 2 µg/ml for WB. |

Telomeric repeat binding factor 2 (TRF2) is a ubiquitously expressed protein that is implicated in the control of telomere length.¹ TRF2, like TRF1, contains a Myb-related DNA binding motif. It binds to duplex TTAGGG repeats and is localized to all human telomeres in metaphase chromosomes.¹ TRF2 is thought to protect chromosome ends by maintaining the correct structure at telomere termini.² The use of mutant forms of TRF2 has implicated a role of TRF2 in the prevention of senescence in primary human cells.² Inhibition of TRF2 results in apoptosis in a subset of mammalian cell types.³

References

1. Broccoli, D., Smogorzewska, A., Chong, L., *et al.* Human telomeres contain two distinct Myb-related proteins, TRF1 and TRF2. *Nat. Genet.* **17**, 231-235 (1997).
2. van Steensel, B., Smogorzewska, A., and de Lange, T. TRF2 protects human telomeres from end-to-end fusions. *Cell* **92**, 401-413 (1998).
3. Karlseder, J., Broccoli, D., Dai, Y., *et al.* p53- and ATM-dependent apoptosis induced by telomeres lacking TRF2. *Science* **283**(5406), 1321-1325 (1999).



Related Products

For a list of related products please visit: www.caymanchem.com/catalog/13791

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, 08/19/2013