

Product Information

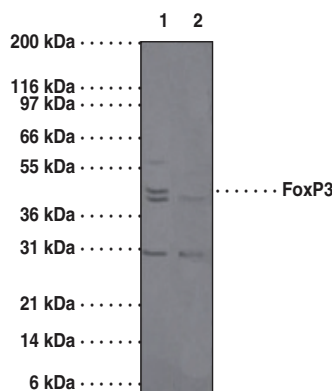


FoxP3/Scurfin Polyclonal Antibody

Catalog No. 13739

- Contents:** This vial contains 100 µg of protein G-purified IgG in 200 µl PBS containing 0.2% gelatin and 0.05% sodium azide.
- Antigen:** Synthetic peptide corresponding to the exon 2 deletion site (i.e. exon 1/exon 3 splice site)
- Host:** Rabbit
- Cross Reactivity:** (+) Bovine, chicken, equine, human, porcine, rat, and Rhesus monkey FoxP3/Scurfin; predicted to react with feline, canine, and murine FoxP3
- Storage:** ≤6 months at 4°C; ≥6 months at -20°C
- Application:** Western blot (WB) and immunohistochemistry. The recommended starting concentration for WB is 2-5 µg/ml.

One of the many immunotolerance mechanisms that the immune system has developed to distinguish between self and non-self antigens are regulatory T-cells or Tregs. Several elegant experiments using transgenic mice and retrovirus-mediated over-expression studies have led to the identification of FoxP3, a transcription factor, as a specific molecular marker essential for the development and function of Tregs. The primary evidence regarding the involvement of FoxP3 in the development of Tregs was provided in patients suffering from IPEX, a rare and fatal human autoimmune disease. The emergence of Tregs and the role of FoxP3 as a critical player in the negative control of various normal and pathological immune responses holds great promise for the development of novel therapies for autoimmune diseases.



Lane 1: FoxP3-transfected cell lysate
Lane 2: Mock-transfected cell lysate

Related Products

FoxP3Δ2 (exon 2 deleted) Specific Monoclonal Antibody (azide free) - Cat. No. 13743 • FoxP3Δ2 (exon 2 deleted) Specific Monoclonal Antibody - Cat. No. 13744

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 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, 02/04/2010