

# PRODUCT INFORMATION



## ROR $\gamma$ Polyclonal Antibody

Item No. 11003

### Overview and Properties

**Contents:** This vial contains 100  $\mu$ g peptide affinity-purified IgG in 200  $\mu$ l PBS, with 0.05% BSA and 0.05% sodium azide.

**Synonyms:** NR1F3, RAR-Related Orphan Receptor C, Retinoid-Related Orphan Receptor  $\gamma$

**Immunogen:** A peptide within the region of amino acids 1-50 of human ROR $\gamma$

**Species Reactivity:** (+) Human and chimpanzee ROR $\gamma$

**Storage:** -20C (as supplied)

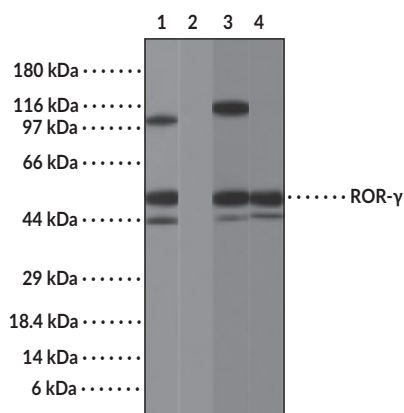
**Stability:**  $\geq$ 6 months

**Host:** Rabbit

**Isotype:** Ig

**Applications:** Immunohistochemistry (IHC) and Western blot (WB); the recommended starting concentration is 10  $\mu$ g/ml for IHC and 0.5-2  $\mu$ g/ml for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

### Image



Lane 1: Jurkat cell lysate (absence of immunizing peptide)  
Lane 2: Jurkat cell lysate (presence of immunizing peptide)  
Lane 3: 3T3 cell lysate  
Lane 4: RAW cell lysate

WARNING  
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA  
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. 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  
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/13/2024

CAYMAN CHEMICAL  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM

# PRODUCT INFORMATION



## Description

---

ROR $\gamma$  is a 560 amino acid novel DNA-binding transcription factor belonging to the ROR/RZR orphan nuclear receptor subfamily. The specific functions of this protein are not known. However, reports suggest that this protein is a ligand activated transcription factor and may be essential for lymphoid organogenesis (lymph nodes and Peyer's patches) and may play an important regulatory role in controlling apoptosis during thymopoiesis. ROR $\gamma$  inhibits the expression of Fas ligand and IL2 and directs the differentiation program of proinflammatory IL-17+ T helper cells. ROR- $\gamma$  shares 50 to 51% sequence identity with ROR $\alpha$  and ROR $\beta$  and is 88% identical to mouse ROR $\gamma$ . It is highly expressed in skeletal muscle and at lower levels in various other tissues.

CAYMAN CHEMICAL  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM