

# Product Information



## ROR $\gamma$ Polyclonal Antibody

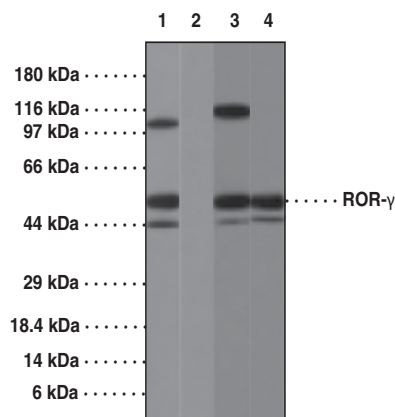
Item No. 11003

<b>Contents:</b>	This vial contains 100 $\mu$ g peptide affinity-purified IgG in 200 $\mu$ l PBS containing 0.05% BSA and 0.05% sodium azide
<b>Synonyms:</b>	NR1F3, RAR-Related Orphan Receptor C, Retinoid-Related Orphan Receptor $\gamma$
<b>Antigen:</b>	A peptide within the region of amino acids 1-50 of human ROR $\gamma$
<b>Host:</b>	Rabbit
<b>Genbank</b>	
<b>Accession No.:</b>	NP_001009598
<b>Isotype:</b>	Rabbit Ig
<b>Cross Reactivity:</b>	(+) Chimpanzee and human ROR $\gamma$ ; predicted to react with bovine, canine, mouse, and rat ROR $\gamma$
<b>Stability:</b>	$\geq$ 6 months at -20°C
<b>Applications:</b>	Immunohistochemistry (IHC) and western blot (WB); the recommended starting concentration for IHC is 10 $\mu$ g/ml and 0.5-2 $\mu$ g/ml for WB.

ROR $\gamma$  is a 560 amino acid 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<sup>+</sup> 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.

### References

1. Kurebayashi, S., Ueda, E., Sakaue, M., *et al.* Retinoid-related orphan receptor  $\gamma$  (ROR $\gamma$ ) is essential for lymphoid organogenesis and controls apoptosis during thymopoiesis. *Proc. Natl. Acad. Sci. USA* **97**(18), 10132-10137 (2000).
2. Ivanov, I.I., McKenzie, B.S., Zhou, L., *et al.* The orphan nuclear receptor ROR $\gamma$ t directs the differentiation program of proinflammatory IL-17<sup>+</sup> T helper cells. *Cell* **126**, 1121-1133 (2006).
3. Eberl, G., Marmon, S., Sunshine, M.-J., *et al.* An essential function for the nuclear receptor ROR $\gamma$ t in the generation of fetal lymphoid tissue inducer cells. *Nature Immuno.* **5**(1), 64-73 (2004).
4. Hirose, T., Smith, R.J., and Jetten, A.M. ROR $\gamma$ : The third member of ROR/RZR orphan receptor subfamily that is highly expressed in skeletal muscle. *Biochem. Biophys. Res. Commun.* **205**(3), 1976-1983 (1994).



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

### Related Products

For a list of related products please visit: [www.caymanchem.com/catalog/11003](http://www.caymanchem.com/catalog/11003)

**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, 05/16/2013

### 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](mailto:custserv@caymanchem.com)

#### Web

[www.caymanchem.com](http://www.caymanchem.com)