

PRODUCT INFORMATION



Transcription Factor Complete PPAR γ Primary Antibody

Item No. 10008897

Overview and Properties

Synonyms: NRIC3, Peroxisome Proliferator-activated Receptor γ
Cross Reactivity: (+) Human PPAR γ
Storage: -20°C (as supplied)
Stability: ≥ 2 years
Host: Rabbit
Applications: PPAR α , δ , γ Complete Transcription Factor Assay Kit

Laboratory Procedures

This antibody has been specifically formulated to be used in PPAR α , δ , γ Complete Transcription Factor Assay Kit (Item No. 10008878). To prepare the primary antibody for use in the PPAR α , δ , γ Complete Transcription Factor Assay Kit, dilute 1:100 in 1X Antibody Binding Buffer (ABB) as outlined below. Add 100 μ l of diluted PPAR γ antibody to each well except the Blank (Blk) wells. Discard any unused diluted primary antibody.

Component	Volume/Well	Volume/Strip	Volume/96-Well Plate
UltraPure Water	89 μ l	712 μ l	8,544 μ l
Transcription Factor Antibody Binding Buffer (10X)	10 μ l	80 μ l	960 μ l
Transcription Factor PPAR γ Primary Antibody	1 μ l	8 μ l	96 μ l
Total Required	100 μ l	800 μ l	9,600 μ l

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 PPAR α , δ , γ Complete Transcription Factor Assay Kit from Cayman (Item No. 10008878). This product may not perform as described if used with other assay reagents or protocols.

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, 10/28/2021

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