

# PRODUCT INFORMATION



## Transcription Factor NF- $\kappa$ B (p50) Primary Antibody

Item No. 10007883

### Overview and Properties

**Synonym:** NF- $\kappa$ B1  
**Host:** Rabbit  
**Cross-reactivity:** (+) Human NF- $\kappa$ B (p50)  
**Storage:**  $\geq 2$  years at  $-20^{\circ}\text{C}$   
**Application:** NF- $\kappa$ B (p50) Transcription Factor Assay Kit

### Description

This antibody has been specifically formulated to be used in NF- $\kappa$ B (p50) Transcription Factor Assay Kit (Item No. 10006912). To prepare primary antibody for use in the NF- $\kappa$ B (p50) Transcription Factor Assay Kit, dilute 1:100 in 1X Antibody Binding Buffer (ABB) as outlined below. Add 100  $\mu\text{l}$  of diluted NF- $\kappa$ B (p50) 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 $\mu\text{l}$	712 $\mu\text{l}$	8,544 $\mu\text{l}$
Transcription Factor Antibody Binding Buffer (10X)	10 $\mu\text{l}$	80 $\mu\text{l}$	960 $\mu\text{l}$
NF- $\kappa$ B (p50) Primary Antibody	1 $\mu\text{l}$	8 $\mu\text{l}$	96 $\mu\text{l}$
Total required	100 $\mu\text{l}$	800 $\mu\text{l}$	9,600 $\mu\text{l}$

### 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 NF- $\kappa$ B (p50) Transcription Factor Assay Kit from Cayman (Item No. 10006912). 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, 06/15/2022

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



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