

Product Information

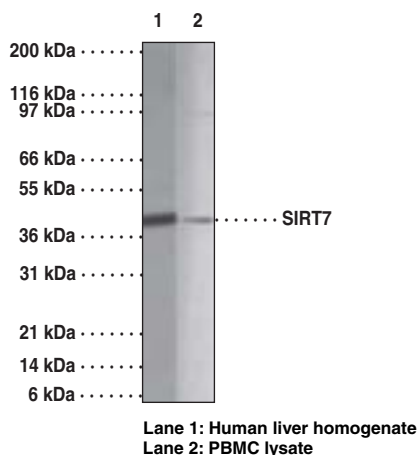


SIRT7 Polyclonal Antibody

Item No. 13477

Contents:	This vial contains 100 µg of protein G-purified IgG in 200 µl PBS containing 0.2% gelatin and 0.05% sodium azide.
Synonyms:	Silent Information Regulator 7, SIR2, Sirtuin 7
Antigen:	Human SIRT7 amino acids 35-51 and 361-377
Host:	Rabbit
Cross Reactivity:	(+) Human SIRT7
Storage:	≤6 months at 4°C; ≥6 months at -20°C
Applications:	The recommended starting dilution for western blot is 1-3 µg/ml.

SIRT7 is a member of a family of proteins called sirtuins (SIR2-like proteins) and are present in prokaryotes and eukaryotes. All SIR2-like proteins have a sirtuin core domain, which contains a series of sequence motifs conserved in organisms ranging from bacteria to humans. Bacterial, yeast, and mammalian sirtuins are able to metabolize NAD⁺ and several act as mono-ADP-ribosyltransferases. The enzymatic function of sirtuins is not yet completely understood but recent reports of histone-activated SIR2-mediated NAD⁺ metabolism and NAD⁺-activated SIR2-mediated histone deacetylation suggest a possible coupled reciprocal activation mechanism involving interactions of SIR2 with NAD⁺ and the N-ε-acetyl-lysine groups of acetylated histone.



Related Products

For a list of related products please visit: www.caymanchem.com/catalog/13477

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, 02/26/2013