

PRODUCT INFORMATION



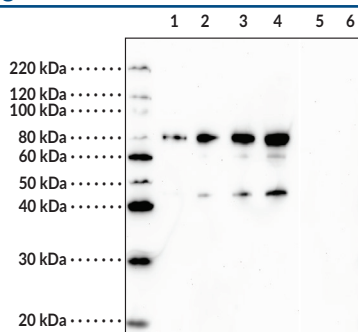
Citrullinated Hsp70 (R155) Polyclonal Antibody

Item No. 24497

Overview and Properties

Contents:	This vial contains 500 µl of peptide affinity purified IgG.
Synonyms:	Heat Shock Protein 70, HspA1A
Immunogen:	Peptide from the internal region of Hsp70 containing a citrulline at residue 155
Cross Reactivity:	(+) Citrullinated Hsp70; (-) Hsp70, other citrullinated Hsp proteins
Species Reactivity:	(+) Human; other species not tested
Uniprot No.:	PODMV8
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥1 year
Storage Buffer:	PBS, pH 7.2, 50% glycerol with 0.02% sodium azide
Host:	Rabbit
Applications:	ELISA and Western blot; the recommended starting dilution is 1:200. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Images



Lane 1: Citrullinated Hsp70 (10 ng)
Lane 2: Citrullinated Hsp70 (25 ng)
Lane 3: Citrullinated Hsp70 (50 ng)
Lane 4: Citrullinated Hsp70 (100 ng)
Lane 5: Hsp70 (200 ng)
Lane 6: Citrullinated Hsp90 (100 ng)

Figure 1. Western blot detection of Hsp proteins using the Citrullinated Hsp70 (R155) Polyclonal Antibody

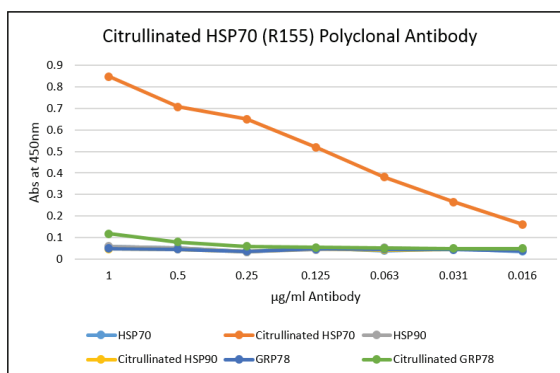


Figure 2. ELISA detection using the Citrullinated Hsp70 (R155) Polyclonal Antibody

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, 03/27/2019

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

Heat shock protein 70s (Hsp70s) are abundant and stress-inducible 70 kDa molecular chaperone proteins encoded by a highly conserved, multigene family.¹ They are monomeric proteins that can be divided into two functional domains: an N-terminal ATPase domain and a substrate binding domain that contains a highly conserved EEVD motif at its C-terminus. Hsp70s are found in the cytosol, nuclei, endoplasmic reticulum, mitochondria, and chloroplasts of eukaryotes, as well as in bacteria. They function as molecular chaperones that assist in a wide range of cellular processes, including refolding of aggregated or misfolded proteins, co- and post-translational folding and assembly of nascent peptides, membrane translocation of secretory and organellar proteins, controlling activity of regulatory nuclear receptors, kinases and transcription factors, as well as cooperativity with the Hsp90 chaperone system in eukaryotes.² The Hsp70 chaperone cycle is ATP-dependent and initiated by transient interaction of the Hsp70 substrate binding domain with hydrophobic regions within a peptide or protein. It consists of an alteration between the low-affinity ATP-bound state with fast rates of substrate exchange and the high-affinity ADP bound state with slow rates of substrate exchange. Hsp70s are subject to a variety of post-translational modifications and their expression is upregulated under conditions of cellular stress and in a variety of disease states. Specifically, citrullinated Hsp70 peptides have been found in the synovial fluid of patients with rheumatoid arthritis.³ Cayman's Citrullinated Hsp70 (R155) Polyclonal Antibody can be used for Western blot and ELISA applications.

References

1. Boorstein, W.R., Ziegelhoffer, T., and Craig, E.A. Molecular evolution of the HSP70 multigene family. *J. Mol. Evol.* **38(1)**, 1-17 (1994).
2. Mayer, M.P. and Bukau, B. Hsp70 chaperones: Cellular functions and molecular mechanism. *Cell Mol. Life Sci.* **62(6)**, 670-684 (2005).
3. Wang, F., Chen, F.F., Gao, W.B., *et al.* Identification of citrullinated peptides in the synovial fluid of patients with rheumatoid arthritis using LC-MALDI-TOF/TOF. *Clin. Rheumatol.* **35(9)**, 2185-2194 (2016).

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