

# PRODUCT INFORMATION



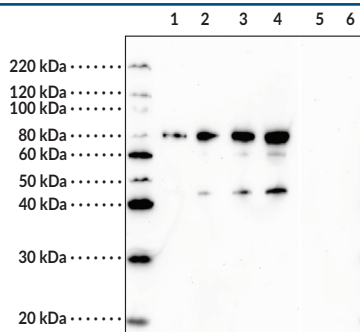
## Citrullinated Hsp70 (R155) Polyclonal Antibody

Item No. 24497

### Overview and Properties

**Contents:** This vial contains 500 µl of peptide affinity purified polyclonal antibody.  
**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.:** P0DMV8  
**Form:** Liquid  
**Storage:** -20°C (as supplied)  
**Stability:** ≥3 years  
**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

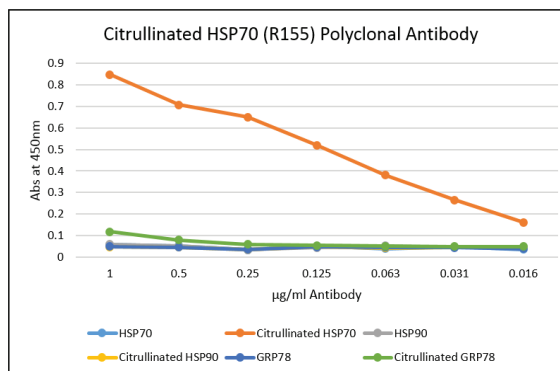


Figure 2. ELISA detection using the Citrullinated Hsp70 (R155) Polyclonal Antibody

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, 11/21/2023

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.<sup>1</sup> 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.<sup>2</sup> 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.<sup>3</sup> 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