

# PRODUCT INFORMATION



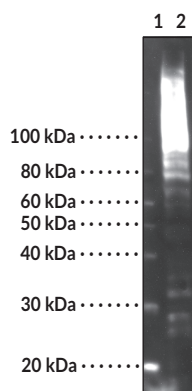
## Multiubiquitin Chain Monoclonal Antibody (Clone FK1)

Item No. 14219

### Overview and Properties

**Contents:** This vial contains 100 µg of affinity-purified monoclonal antibody.  
**Immunogen:** Crude preparation of polyubiquitinated-lysozyme  
**Cross Reactivity:** (+) Polyubiquitinated conjugates; (-) Free ubiquitin and monoubiquitinated conjugates  
**Species Reactivity:** (+) Species independent  
**Form:** Liquid  
**Storage:** -20°C (as supplied)  
**Stability:** ≥1 year  
**Storage Buffer:** 10 mM phosphate buffer, pH 7.4, with 0.15 M sodium chloride, and 0.1% sodium azide  
**Clone:** FK1  
**Host:** Mouse  
**Isotype:** IgM  
**Applications:** Western blot (WB); the recommended starting dilution is 1:300. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.  
**Special Conditions:** Milk based blocking reagents are not recommend. 1% BSA in PBS or TBS polysorbate 20 is recommended as a blocking buffer.

### Image



Lane 1: MW Markers  
Lane 2: HeLa cell extract (20 µg)

WB analysis using Multiubiquitin Chain Monoclonal Antibody (Clone FK1) at a dilution of 1:300.

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, 01/21/2021

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

# PRODUCT INFORMATION



## Description

---

Ubiquitin is a small protein composed of 76 amino acids and having a molecular weight of 8.6 kDa. It exists in all eukaryotic cells and is responsible for numerous functions in the cell by post-translational modifications of substrate proteins (ubiquitination). This process is important to numerous cellular processes including, but not limited to, apoptosis, cell cycle and division, DNA transcription, immune response, modulation of cell surface receptors, and viral infections. Clone FK1 reacts with polyubiquitinated conjugates, but does not react with free ubiquitin or monoubiquitinated conjugates.<sup>1,2</sup> Cayman's Multiubiquitin Chain Monoclonal Antibody (Clone FK1) can be used for Western blotting applications in species independent samples.

## References

---

1. Takada, K., Nasu, H., Hibi, N., *et al.* Serum concentrations of free ubiquitin and multiubiquitin chains. *Clin. Chem.* **43(7)**, 1188-1195 (1997).
2. Fujimuro, M., Sawada, H., and Yokosawa, H. Production and characterization of monoclonal antibodies specific to multi-ubiquitin chains of polyubiquitinated proteins. *FEBS Lett.* **349(2)**, 173-180 (1994).

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