

# PRODUCT INFORMATION



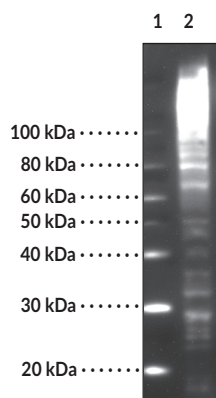
## Multiubiquitin Chain Monoclonal Antibody (Clone FK2)

Item No. 14220

### Overview and Properties

<b>Contents:</b>	This vial contains affinity purified monoclonal antibody.
<b>Immunogen:</b>	Crude preparation of polyubiquitinated-lysozyme
<b>Cross Reactivity:</b>	(+) Mono- and polyubiquitinated conjugates; (-) free ubiquitin
<b>Form:</b>	Liquid
<b>Storage:</b>	-20°C (as supplied)
<b>Stability:</b>	≥1 year
<b>Storage Buffer:</b>	10 mM phosphate buffer, pH 7.4, with 0.15 M sodium chloride and 0.1% sodium azide
<b>Concentration:</b>	1 mg/ml (100 µg vial) or 10 mg/ml (500 µg vial)
<b>Clone:</b>	FK2
<b>Host:</b>	Mouse
<b>Isotype:</b>	IgG <sub>k1</sub>
<b>Applications:</b>	ELISA, immunohistochemistry, and western blot (WB); the recommended starting dilution for WB is 1:100-1:1,000. Other applications were not attempted and therefore optimal working dilutions should be determined empirically.

### Image



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

WB analysis using Multiubiquitin Chain Monoclonal Antibody (Clone FK2) at a 1:1,000 dilution.

**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, 01/20/2021

**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 kD. 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 FK2 is the most widely used ubiquitin antibody. This clone reacts with mono- and polyubiquitinated conjugates, but does not react with free ubiquitin.<sup>1,2</sup>

## 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