

PRODUCT INFORMATION



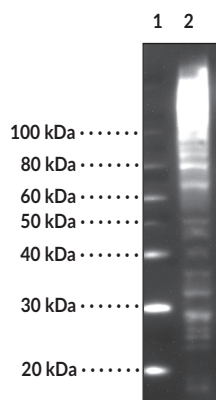
Multiubiquitin Chain Monoclonal Antibody (Clone FK2)

Item No. 14220

Overview and Properties

Contents:	This vial contains affinity-purified monoclonal antibody.
Immunogen:	Crude preparation of polyubiquitinated-lysozyme
Cross Reactivity:	(+) Mono- and polyubiquitinated conjugates; (-) free ubiquitin
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥3 years
Storage Buffer:	PBS, pH 7.2, with 0.09% sodium azide
Concentration:	1 mg/ml (100 µg vial) or 10 mg/ml (500 µg vial)
Clone:	FK2
Host:	Mouse
Isotype:	IgG _{k1}
Applications:	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, 11/14/2023

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

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.^{1,2}

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