

[pThr¹³⁴²]DNA-topoisomerase II α monoclonal antibody (3D4)

DNA-topoisomerase II, also known as Topo II or Top 2, is a multifunctional nuclear enzyme required during DNA replication, chromosome disjunction at mitosis, and other DNA-related activities because of its ability to alter DNA supercoiling. In mammalian cells, Topo II consists of two isozymes, Topo II-alpha (170kDa) and Topo II-beta (180kDa) and expression and localization of each isoform are distinct and stage specific during the cell cycle. Topo II-beta is expressed constitutively throughout the cell cycle, whereas the expression of Topo II-alpha is cell cycle-regulated, peaking in G2 to M phase and declining to a minimal level at the end of M phase. DNA-topoisomerase II is the target of some of the most active drugs used in chemotherapy and has been shown to have a critical role in neural development.

This antibody is covered by our [Worry-Free Guarantee](#).

Ordering Information

[Order Online »](#)

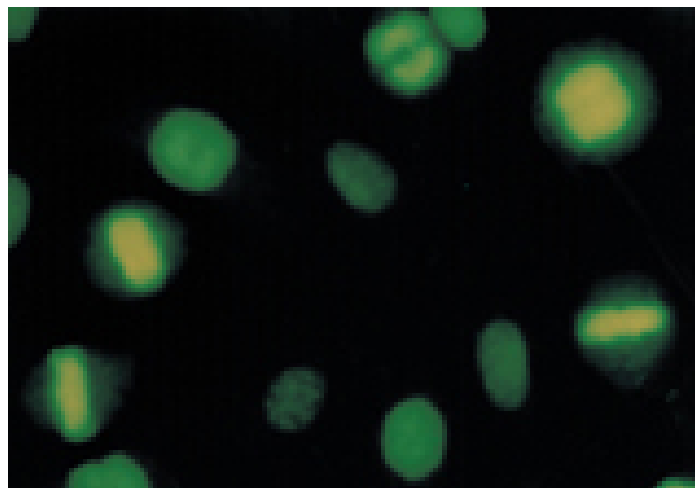
| | |
|-----------------|-------|
| ADI-KAM-CC220-E | 100µg |
|-----------------|-------|

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of Jurkat cells using DNA-Topoisomerase IIα (pThr1342) mAb (3D4).



Immunocytochemistry analysis of human HEP-2 cells stained with DNA-Topoisomerase IIα (pThr1342) mAb (3D4).

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application ICC, WB

Application Notes Detects a band of ~170kDa by Western blot.

Clone 3D4

Formulation Liquid. In PBS, pH 7.2, containing 50% glycerol.

Host Mouse

Immunogen Synthetic peptide corresponding to a portion of human Topoisomerase II α phosphorylated at Tyr1342.

Isotype IgG1

Purity Detail Protein A affinity purified.

Recommendation Immunocytochemistry (10 μ g/ml)Western Blot (1-10 μ g/ml, colorimetric)Suggested
Dilutions/Conditions dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Species Reactivity Human

UniProt ID P11388 (α isozyme), Q02880 (β isozyme)

Worry-free Guarantee This antibody is covered by our [Worry-Free Guarantee](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)