

Tyrosyl-DNA Phosphodiesterase 1 (human) (recombinant) (His)

TDP1 is catalytically active and will remove a terminal 3' phosphotyrosyl group from a specifically designed oligo target site (oligo-Tyr). Additionally, TopoGEN scientists have demonstrated that TDP1 will release topo I covalently attached to genomic DNA isolated from cells treated with camptothecin using TopoGEN's *In Vivo* Link Kit ([Prod. No. TOP-1021](#)).

Ordering Information

[Order Online »](#)

TOP-TG2004H-1	250U
---------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	Avoid freeze/thaw cycles.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name TDP1

Specific Activity One unit of TDP1 is defined as the amount of enzyme which can release the terminal Tyr from 20ng of substrate (oligo-Tyr) in 30 min. at 37°C.



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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