

PLK1 (human), (recombinant) (His- tag)

PLK localizes to the mitotic spindle and microinjection of PLK into quiescent NIH 3T3 cells induces them to enter mitosis. PLK1 is elevated in many types of cancer and when depleted by siRNA, cell proliferation is inhibited and viability decreased.

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

| | |
|----------------|-----|
| BML-SE466-0005 | 5µg |
|----------------|-----|

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Polo-like kinase 1, STPK13, Serine/threonine protein kinase 13

Application Notes Useful for kinetic and functional studies, phosphorylation of target substrates, drug screening.

MW 70 kDa

Source Produced in insect cells. Full length active PLK1 is fused at the N-terminus to a His-tag. Produced in a baculovirus expression system.

UniProt ID P53350

Last modified: May 29, 2024



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)