

Flt1 (human), (recombinant) (His- tag)

Flt1 (FMS-related Tyrosine Kinase 1, VEGFR1) is implicated in preeclampsia, control of metastasis, and in wound repair.

Ordering Information

[Order Online »](#)

BML-SE445-0010	10µg
-----------------------	------

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

FMS-related tyrosine kinase 1, VEGFR-1, Vascular endothelial growth factor receptor 1

Application Notes

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

MW

68 kDa

Source

Produced in insect cells. Active Flt1 (aa 783- end).
Produced in a baculovirus expression system.

UniProt ID

Q8QHL3

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)