

CD105 (human), mAb (SN6/N1-3A1)

Ordering Information

[Order Online »](#)

ANC-326-030	Biotin - 100µg
ANC-326-020	Purified - 100µg
ANC-326-820	Preservative Free - 100µg
ANC-326-040	FITC - 120 tests
ANC-326-050	R-PE - 120 tests
ANC-326-060	APC - 120 tests

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Endoglin

Application Flow Cytometry, IHC

Application Notes Functional Application (augments binding of TGF β 1 to CD105 expressing leukemia cells).

Clone SN6/N1-3A1

Host Mouse

Immunogen Membrane preparation of human B-lineage leukemia cells.

Isotype IgG1

Species Reactivity Horse, Human, Monkey, Porcine

Specificity Recognizes human CD105 (endoglin) as a homodimer.

Technical Info / Product Notes For the Original Manufacturer's Data Sheet please click on the corresponding product number:

[ANC-326-020](#)

[ANC-326-030](#)

[ANC-326-040](#)

[ANC-326-050](#)

[ANC-326-060](#)

[ANC-326-820](#)



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-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-de@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