

Annexin V (human), (recombinant) (FITC conjugate)

FITC-labelled recombinant human annexin V fusion protein (produced in *E. coli*) shows bright green fluorescence (Ex(max): 488nm; Em(max): 530nm).

Citations: 4

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-209-256-T100	100 tests
ALX-209-256-3100	3x100 tests

Manuals, SDS & CofA

[View Online »](#)

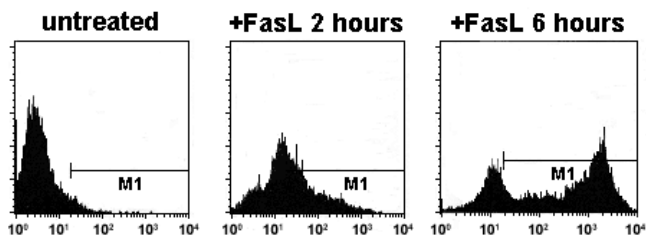
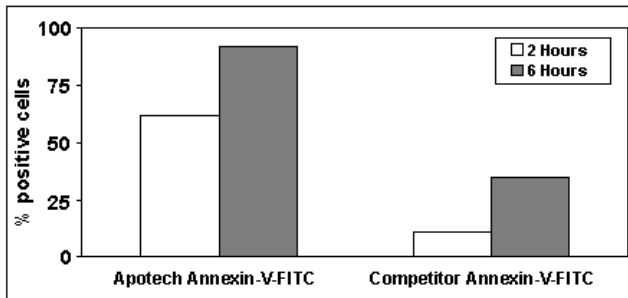
A**B**

Figure A: FACS histogram for the fluorescence intensity measured in FL-1 (FITC) channel. Jurkat cells were incubated for 2 or 6 hours with medium (untreated) or with 100ng/ml rhsFasL in the presence of 1 μ g/ml Enhancer for Ligands (rhsFasL Set, Prod. No. [ALX-850-014](#)).

Figure B: Comparison of ALEXIS and Competitor Annexin-V. Jurkat cells were incubated for 2 or 6 hours with medium (background) or with 100ng/ml rhs FasL in the presence of 1 μ g/ml Enhancer for Ligands (rhs FasL Set, Prod. No. 850-014). Cells were stained with Annexin-V-FITC (Prod. No. 209-256 and competitor”

Handling & Storage

Use/Stability	Stable for at least one year after receipt when stored at +4°C.
Handling	Protect from light. Do not freeze.
Long Term Storage	+4°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application Notes	Detection of apoptotic cells by flow cytometry.
Concentration	1mg/ml
Formulation	Liquid. In PBS, pH 7.4, containing 0.02% sodium azide.
MW	~52kDa (SDS-PAGE).
Purity	≥90%
Purity Detail	Affinity purified.
Quantity	Sufficient for 100 or 300 tests. 1 test is equivalent to 1µg of the conjugated protein in an experimental volume of 0.4 ml.
Source	Produced in <i>E.coli</i> .
Specificity	Binds to phosphatidylserine (PS).
UniProt ID	P08758

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-
fr@enzolifesciences.com](mailto:info-fr@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)