

Annexin V (human), (recombinant) (R-PE conjugate)

R-PE-labelled recombinant human annexin V is suitable for dual staining experiments to detect phosphatidylserine whilst cell-surface protein can be detected with a suitable MAb (FITC-labelled or labelled secondary antibody). The fluorochrome R-PE (phycoerythrin) shows red fluorescence (Ex(max): 495-564nm; Em(max): 576nm). Exhibits anti-phospholipase activity.

Ordering Information

[Order Online »](#)

ALX-209-2510	100 tests
--------------	-----------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

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.
Formulation	Liquid. In 50mM TRIS, pH 7.4, containing 100mM sodium chloride, 1% BSA and 0.02% sodium azide.
MW	~35.8kDa.
Purity	≥89% (SDS-PAGE, HPLC)
Quantity	20 tests/100µl; 100 tests/500µl.
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

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