

Alkaline Phosphatase Staining Kit

For visualization of ALP activity in cultured cells under a microscope.

Alkaline Phosphatase (ALP) is an enzyme found in various tissues throughout the body, including the liver, bone, kidney, and intestine. It plays a crucial role in catalyzing the hydrolysis of phosphate groups from various molecules including nucleotides and proteins. ALP is essential for many physiological processes, including bone mineralization, nucleic acid metabolism, and cellular signaling. AkrivisBio's Alkaline Phosphatase utilizes the BCIP/NBT protocol based on indoxyl substrates. In the stain, BCIP is hydrolyzed to an indoxyl, which reduces the NBT to a darkly colored formazan.

Ordering Information

[Order Online »](#)

AKR-COS-0105	1Kit
--------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Species Reactivity Species independent

Technical Info / Product Notes For the Original Manufacturer's data sheet please [click here](#).



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