

Beta-galactosidase Staining Kit

For measuring β -Galactosidase in various samples.

The E. coli LacZ gene is a commonly used reporter gene for testing the expression efficiency of vector mediated gene transfer. It is also used frequently to study gene promoter regulation. LacZ encodes β -galactosidase, an almost perfectly ideal choice for expression due to being very stable, resistant to proteolytic degradation, able to utilize multiple substrates and easily assayed in situ. AkrivisBio's β -Galactosidase Staining Kit is a simple, sensitive assay, which utilizes X-Gal a colorless substrate, resulting in an intense blue insoluble indigo product easily visualized under microscopy.

Ordering Information

[Order Online »](#)

AKR-COS-0104	1Kit
--------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°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