

Phenylalanine Assay

A plate based fluorometric assay to measure phenylalanine in various sample types with a sensitivity in the mid to low picomole range.

Phenylalanine is one of the essential amino acids that we are unable to synthesize. Beyond its obvious significance as an important building block for protein synthesis, phenylalanine is also a precursor dopamine, norepinephrine, and epinephrine as well as a precursor for tyrosine which itself is a precursor to thyroid hormones. PKU is a genetic disorder characterized by the inability to degrade phenylalanine resulting in its accumulation causing neurological and cognitive impairment, at least partly due to competing for tryptophan transport channels resulting in decreased serotonin. AkrivisBio's Phenylalanine Assay is a simple, sensitive means of measuring phenylalanine in a variety of biological samples with a sensitivity in the mid to low picomole range.

Ordering Information

[Order Online »](#)

AKR-MA-0147	100 wells
-------------	-----------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Fluorescent detection

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@enzolifesciences.com
usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info@enzolifesciences.com
eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info@enzolifesciences.com
be@enzolifesciences.com

France
Phone: +33 472 440 655
info@enzolifesciences.com
fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info@enzolifesciences.com
de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info@enzolifesciences.com
uk@enzolifesciences.com