

# Iron Assay

A plate-based colorimetric assay to measure 0-10 nanomoles of iron levels in samples.

Iron is an essential element involved in many physiological processes including erythropoiesis, immune function, mitochondrial function, electron transport, DNA synthesis and enzymatic reactions, which require iron as a cofactor. The most common structures in which iron is found are the iron-sulfur centers and various hemes. Unlike most other ions whose levels are regulated by both absorption and excretion, iron levels in vivo are controlled solely by absorptive processes. AkrivisBio's Iron Assay is a simple and direct method for measuring (Fe<sup>2+</sup>) and/or ferric (Fe<sup>3+</sup>) iron in samples. In this assay, iron compounds present in samples are optionally reduced to ferrous form (Fe<sup>2+</sup>), which then reacts with Ferene S to produce a stable chromogen measured at 593 nm. The assay can measure iron in the range of 0-10 nanomoles in a variety of samples.

## Ordering Information

[Order Online »](#)

AKR-MA-0103

100 wells

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Application Colorimetric 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-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-be@enzolifesciences.com](mailto:info-be@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)