

Branched Chain Amino Acid Assay

A plate based colorimetric assay to measure BCAA's in a variety of sample types with a sensitivity down into the low nanomole range vs normal serum concentrations in the

50-300 nmole/ml range. The branched-chain amino acids, comprising leucine, isoleucine, and valine, are deemed essential, and must be obtained through diet as the body cannot produce them independently. BCAAs are particularly important for muscle health and growth. Leucine, in particular, acts as a powerful stimulator of protein synthesis and insulin secretion. BCAAs also act as an energy source during periods of intense exercise, preserving muscle mass and preventing their breakdown. BCAAs are unique in their ability to bypass the liver and be directly utilized by muscles for energy production. An imbalance in BCAA metabolism is linked to various health conditions like insulin resistance, obesity, and muscle wasting. Some of this appears to occur through mTORC impairment. On the other hand, maintaining an appropriate BCAA ratio is associated with improved insulin sensitivity and a reduced risk of chronic diseases. AkrisBio's BCAA Assay is a simple direct method of measuring the BCAA's in a variety of sample types with a sensitivity down into the low nanomole range vs normal serum concentrations in the 50-300 nmole/ml range.

Ordering Information

[Order Online »](#)

AKR-MA-0157	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 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

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