

β -Alanine

β -Alanine is a common non-essential beta amino acid formed by degradation of dihydrouracil and carnosine. Levels of β -alanine limit the amount of carnosine, and it has been used frequently as a workout supplement to increase carnosine concentrations in muscle. This assists in building lean muscle mass, and can also improve physical functioning and performance.

Materials Provided

Catalog Number:	92-1227
Quantity Supplied:	1 vial(s), 10 mg per vial

Description

DiscoverX control ligands have been functionally tested and validated for optimal use with all cell line targets and Assay Ready kits.

Product Information

Molecular Weight:	89.09 g/mol
Source:	Synthetic
Purity:	$\geq 98\%$ TLC
Endotoxin Level:	N/A
Formulation:	N/A
Storage:	RT Please avoid multiple freeze/thaw cycles.

Reagent Preparation

To avoid condensation, equilibrate the vial to ambient temperature before opening.

Stock Concentration:	100 mM
Reconstitution Volume:	1.122 mL
Reconstitution Solvent:	H ₂ O

For Research Use Only

Ordering: +1.510.979.1415 option 4 or e-mail CustomerServiceDRX@eurofins.com
Technical support: +1.510.979.1415 option 5 or e-mail DRX_SupportUS@eurofinsUS.com
General product information: www.discoverx.com