

FICZ

FICZ (5,11-dihydro-indolo[3,2-b]carbazole-6-carboxaldehyde, 6-Formylindolo[3,2-b]carbazole) is a high affinity aryl hydrocarbon receptor (AhR) agonist that induces transient expression of cytochrome P450-1A1 (CYP1A1) in vitro and plays roles in immune response, control of inflammation, cell-cycle regulation, and development and maturation of many tissues.

Materials Provided

Catalog Number:	92-1329
Quantity Supplied:	1 vial(s), 1 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:	284.31 g/mol
Source:	Synthetic
Purity:	>95%
Endotoxin Level:	N/A
Formulation:	N/A
Storage:	-20°C Please avoid multiple freeze/thaw cycles.

Reagent Preparation

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

Stock Concentration:	10 mM
Reconstitution Volume:	0.352 ml
Reconstitution Solvent:	DMSO

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