

6-Formylindolo(3,2-b)carbazole

Ah receptor ligand

A tryptophan-derived, high affinity aryl hydrocarbon receptor ligand ($K_d = 7 \times 10^{-11}$ M). Proposed to be an endogenous AhR ligand. Induces rapid and transient expression of cytochrome P450-1A1 (CYP1A1) in human cells at 100 pM.

Citations: 51

[View Online »](#)

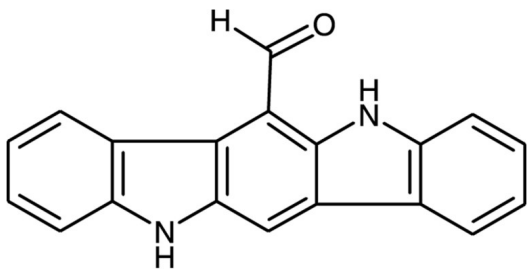
Ordering Information

[Order Online »](#)

BML-GR206-0100	100µg
----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store solutions at -20°C for up to one month.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	FICZ
Appearance	Yellow solid.
CAS	172922-91-7
Couple Target	Ah receptor
Couple Type	Ligand
Formula	$C_{19}H_{12}N_2O$
MW	284.3
MeltingPoint	>300°C
Purity	≥95% (HPLC)
Solubility	Soluble in DMSO (10mg/ml).

Last modified: May 29, 2024



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