

MDBN

3,4-Methylenedioxy- β -nitrostyrene

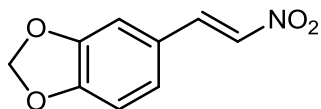
Cat. # SI9770

Background: An irreversible inhibitor of p97/vasolin ($IC_{50}=1.7 \mu M$) identified in a cell-based protein degradation assay. Cell permeable.

Application: Cell-based or in vitro characterization of p97/vasolin function

Product Information:

CAS No.	1485-00-3
Purity:	>98% (HPLC); NMR (Conforms)
Molecular Weight:	193.2
Physical State:	Yellow Powder
Quantity:	25 mg
Solubility:	May be dissolved in DMSO (50 mg/ml)
Storage:	Store desiccated as supplied at ambient temperature for up to 2 years. Store solutions at -20°C for up to 1 month.



Formula: $C_9H_7NO_4$

References

1. Chou T. F. and Deshaies R. *et al.* Quantitative cell-based protein degradation assays to identify and classify drugs that target the ubiquitin-proteasome system. *J. Biol. Chem.* 2011 **286**:16546

All products are for research use only • Not intended for human or animal diagnostic or therapeutic uses
Copyright © 2011 LifeSensors, Inc. All Rights Reserved