

PRODUCT: PathwayReady™ MAP Kinase Signaling Inhibitor Panel

CATALOG #: K862-7

AMOUNT: 1 set

STORAGE: Store at -20°C. Protect from air and light

DESCRIPTION: A convenient set consisting of seven inhibitors useful for the study of MAP Kinase signaling pathway. The seven inhibitors are shown in the following table:

Product	Cat. No.	Size	Biological Activity
PD 184352 (CI-1040)	1585-1	1 mg	MEK1/2 inhibitor
PD 98059	1666-5	5 mg	MEK1 inhibitor
SB 203580	1712-1	1 mg	p38 MAP Kinase inhibitor
Sorafenib	1594-25	25 mg	Raf, PDGF, VEGF inhibitor
SP 600125	1669-5	5 mg	JNK inhibitor
U0126	1668-5	5 mg	MEK1/2 inhibitor
ZM 336372	1789-1	1 mg	c-Raf inhibitor

HANDLING: Do not take internally. Wear gloves and mask when handling the product! Avoid contact by all modes of exposure.

USAGE: *FOR RESEARCH USE ONLY! Not to be used in humans.*

RELATED PRODUCTS:

- PD 184352 (CI-1040) (Cat. No. 1585-5, 25)
- PathwayReady™ PI3-K/Akt/mTOR Signaling Pathway Inhibitor Panel (Cat. No. K857-12)
- PD 98059 (Cat. No. 1666-5)
- EZSolution™ PD 98059 (Cat. No. 1783-5)
- PD-0325901 (Cat. No. 1643-2)
- SB 203580 (Cat. No. 1712-1, 5)
- EZSolution™ SB 203580 (Cat. No. 1786-1)
- U0126 (Cat. No. 1668-5)
- Sorafenib (Cat. No. 1594-25, 100)
- VX-702 (Cat. No. 1598-1,5)

