

PRODUCT: DiscoveryPak™ PI 3-Kinase Inhibitor Set

CATALOG #: K856-5

AMOUNT: 1 set

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

DESCRIPTION: A convenient set consisting of five phosphoinositide 3-kinase (PI 3-kinase) inhibitors with diverse biological activities. The five products are shown in the following table:

Product	Cat. No.	Size	Biological Activity
IC87114	1661-1	1 mg	PI3-K δ inhibitor
LY 294002	1667-5	5 mg	Inhibits PI3-Ks
PI-103	1728-1	1 mg	PI3-K α inhibitor
PI3-K γ Inhibitor, AS-605240	1780-1	1 mg	PI3-K γ Inhibitor
TGX-221	1781-1	1 mg	PI3-K β 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:

- BEZ235 (NVP-BE235) (Cat. No. 1626-5, 25)
- GDC-0941 bismesylate (Cat. No. 1623-1, 5)
- Deforolimus (Cat. No. 1587-5, 25)
- IC87114 (Cat. No. 1661-1)
- LY 294002 (Cat. No. 1667-5, 25)
- mTOR Inhibitor, Ku-0063794 (Cat. No. 1779-1,5)
- PI-103 (Cat. No. 1728-1,5)
- PI3-K γ Inhibitor, AS-605240 (Cat. No. 1780-1,5)
- PP242 (Cat. No. 1658-1)
- Rapamycin (Cat. No. 1568-5, 50)
- Temsirolimus (Cat. No. 1600-5, 25)
- TGX-115 (Cat. No. 1660-1)
- TGX-221 (Cat. No. 1781-1,5)
- Wortmannin (Cat. No. 1670-1)

