

**PRODUCT: PathwayReady™ JAK/STAT Signaling Inhibitor Panel**

**CATALOG #:** K864 -8

**AMOUNT:** 1 set

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

**DESCRIPTION:** A convenient set consisting of eight inhibitors useful for studying JAK/STAT signaling pathway. The 8 inhibitors are shown in the following table:

Product	Cat. No.	Size	Inhibits
AG 490	1570-5	5 mg	JAK2 and JAK3
CP-690550	1622-5	5 mg	JAK3
Lestaurtinib	1805-500	500 µg	JAK2/STAT5
LLL12	1792-5	5 mg	STAT3
Pimozide	1887-5	5 mg	STAT5
Stat3 Inhibitor, S3I-201	1775-5	5 mg	STAT3
Stat3 Peptide Inhibitor, Cell-Permeable	1778-1	1 mg	STAT3
WP1066	1809-5	5 mg	JAK2, STAT3

**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:**

- DiscoveryPak™ EGFR Tyrosine Kinase Inhibitor Set (Cat. No. K858-5)
- DiscoveryPak™ Receptor Tyrosine Kinase (RTK) Inhibitor Set (Cat. No. K859-8)
- DiscoveryPak™ PI3- Kinase Inhibitor Set (Cat. No. K856-5)
- PathwayReady™ EGFR Signaling Inhibitor Panel (Cat. No. K863-14)
- PathwayReady™ MAP Kinase Signaling Inhibitor Panel (Cat. No. K862-7)
- PathwayReady™ PI3-K/Akt/mTOR Signaling Inhibitor Panel (Cat. No. K857-12)

