

## DiscoveryPak™ Hedgehog Signaling Pathway Inhibitors Set

**CATALOG #:** K868-6

**AMOUNT:** 1 set

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

**DESCRIPTION:** A convenient set consisting of five Hedgehog (Hh) signaling pathway inhibitors and a negative control. The six products are shown in the following table:

Product	Cat. No.	Size	Biological Activity
Cyclopamine-KAAD	1910-50	50 µg	Hh pathway inhibitor
GANT58	1812-5	5 mg	Hh pathway inhibitor/Gli antagonist
GDC-0449	1890-5	5 mg	Hh pathway inhibitor
Hh Signaling Pathway Antagonist	1659-1	1 mg	Hh pathway inhibitor
JK 184	1726-1	1 mg	Hh pathway inhibitor
Tomatidine hydrochloride	1893-25	25 mg	Negative control

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

### RELATED PRODUCTS:

- Cyclopamine (Cat. No. 1578-5)
- Cyclopamine-KAAD (Cat. No. 1910-50)
- GANT58 (Cat. No. 1812-5, 25)
- GANT61 (Cat. No. 1892-5)
- GDC-0449 (Cat. No. 1890-5, 25)
- Hh Signaling Pathway Antagonist (Cat. No. 1659-1)
- JK-184 (Cat. No. 1726-1)
- Purmorphamine (Cat. No. 1672-5)
- Tomatidine hydrochloride (Cat. No. 1893-25)

**FOR RESEARCH USE ONLY! Not to be used on humans.**