

PRODUCT: DiscoveryPak™ DPP-IV Inhibitors Set**CATALOG #:** K890-5**AMOUNT:** 5 inhibitors**STORAGE:** Store at -20°C. Protect from light.**DESCRIPTION:** A convenient set of five (5) small molecule Dipeptidyl IV (DPP-IV) inhibitors (as shown in the table below).

Product	Cat. No.	Size
Linagliptin	2240	50 mg
Saxagliptin	9528	50 mg
Sitagliptin phosphate Monohydrate	1757	100 mg
Trelagliptin Succinate	2798	5 mg
Vildagliptin	2188	10 mg

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

Dipeptidylpeptidase IV, Human Plasma (**Cat. No. 4709-10**)
Dipeptidylpeptidase IV, Human Recombinant (**Cat. No. 4710-10, 50, 1000**)
Dipeptidylpeptidase IV Inhibitor, K 579 (**Cat. No. 1963-1, 5**)
Dipeptidylpeptidase IV Inhibitor, NVP DPP 728 (**Cat. No. 1964-1, 5**)
Diprotin A (**Cat. No. 2191-5, 25**)
Diprotin B (**Cat. No. 2192-5, 25**)
DPP4 Activity Assay Kit (**Cat. No. K779-100**)
DPP4 Inhibitor Screening Kit (**Cat. No. K780-100**)
Linagliptin (**Cat. No. 2240-50, 250**)
Sitagliptin Phosphate Monohydrate (**Cat. No. 1757-100, 1G**)
Vildagliptin (**Cat. No. 2188-10, 50**)

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