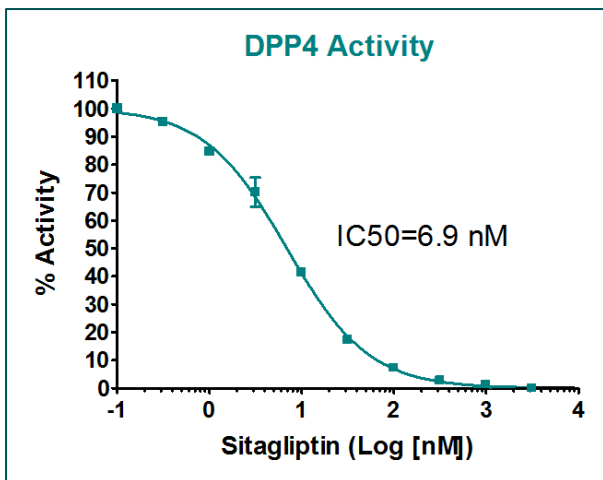
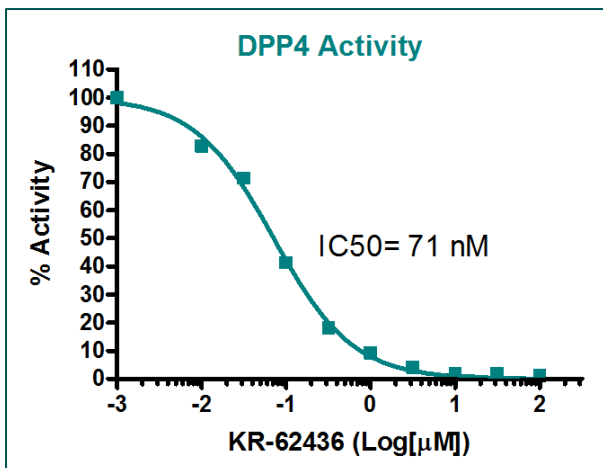


DPP and other Peptidase Profiling Services



Sample IC₅₀ assays for DPP4 activity, using KR-62436 (top image) or sitagliptin (bottom)

DPP (Dipeptidyl Peptidase) Screening & Profiling:

- DPP3
- DPP4
- DPP7
- DPP8
- DPP9
- FAP
- POP

Other Peptidase Screening & Profiling:

- Cathepsin B
- Cathepsin D
- Cathepsin E
- Cathepsin F
- Cathepsin L
- Cathepsin S
- Cathepsin V

Other Screening/Profiling Services:

HDACs (all 11 available)

Histone Demethylases (4 available)

Histone Methyltransferases (14 available)

SIRTs (all 7 available)

PARPs (8 available)

PDEs (23 available)

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082**, Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com