

Data Sheet

DPP8

Human, recombinant, N-terminal GST tag

Catalog #: 80080

Lot#: 120827-GC **Conc.:** 1.12 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
 110 mM NaCl, 2.2 mM KCl, and 20%
 glycerol.

Stability: At least 6 months at -80°C

References:

1. Lee, H.J. *et al.*, *J. Biol. Chem.* **281** (50), 38653-38662 (2006).
2. Yu, D.M. *et al.*, *FEBS J.* **273** (11), 2447-2460 (2006).

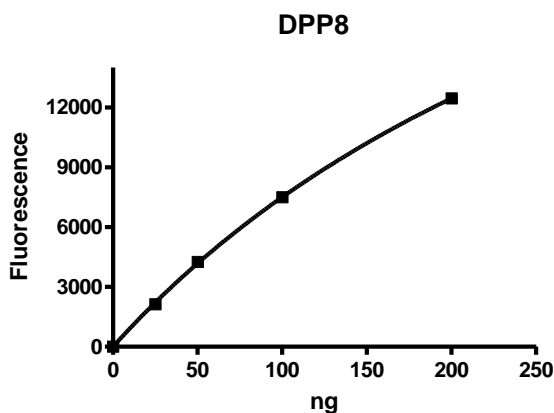
Description: Human DPP8 also known as, Dipeptidyl-peptidase 8, GenBank Accession No. NM_197960, a.a. full length with N-terminal GST tag, expressed in a Baculovirus infected Sf9 cell expression system, MW=129.3 kDa.

Specific Activity: ≥ 560 pmol/ min/ μg
Assay Conditions: Ala-Pro-AMC was used as a substrate. Reaction was done in a buffer containing 10 mM Tris (pH 7.4), 10 mM MgCl_2 , 0.05 % Tween-20 and 10 μM substrate for 30 minutes at room temperature.

Application: Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance

Specific Activity:



4-20% SDS-PAGE
Coomassie staining

Lane 1:
 2 μg DPP8
Lane 2:
 Protein Marker

MW: 129.3 kDa.
Purity: $\geq 80\%$

