

## FLT3 (D835Y)

(FL cytokine receptor [Precursor], Tyrosine-protein kinase receptor FLT3, Stem cell tyrosine kinase 1, STK-1, Flk-2)

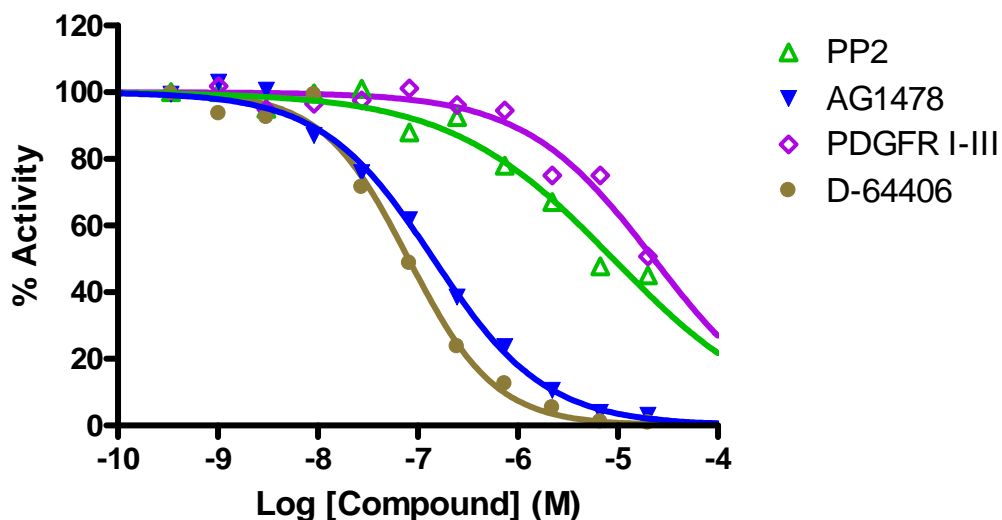
- **CAT#:** FLT3 (D835Y)
- **Enzyme:** Human FLT3 (D835Y)
- **Substrate:** Peptide substrate, [EAIYAAPFAKKK], 5  $\mu$ M  
ATP 10  $\mu$ M
- **Reaction:**  

$$\text{Substrate} + [\gamma\text{-}^{33}\text{P}]\text{-ATP} \xrightarrow{\text{Enzyme}} \text{}^{33}\text{P-Substrate} + \text{ADP}$$
- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	< 0.1
PP2	9,000
AG1478	143
PDGFR I-III	22,800
D-64406	80.8

- **Scientific Information:** P36888 (<http://www.expasy.org/uniprot/P36888>)

### Reference compound IC50 for FLT3 (D835Y)



Service performed  
using kinases  
under license from  **invitrogen**<sup>™</sup>  
part of *life* technologies<sup>™</sup>