

Kinase mutant panel performed with ³³PanQinase™ assay technology.

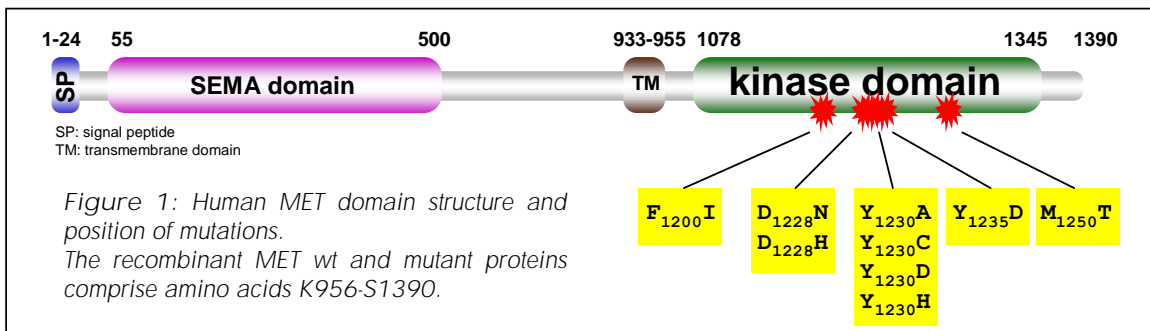
➤ MET Kinase _{K956-S1390} Mutant Panel

met proto-oncogene

Synonyms: c-MET, HGFR

MET plays a pivotal role in tumor growth, metastasis and angiogenesis. Many pathophysiological and oncogenic MET mutants have been described. Furthermore, several MET mutations confer resistance against therapeutical MET kinase inhibitors^{1,2}.

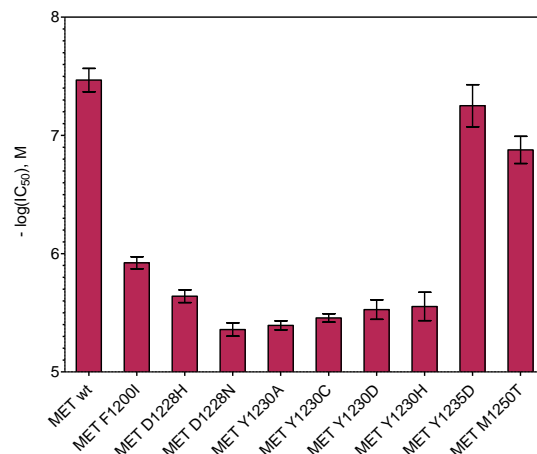
MET wildtype (wt) and nine pathophysiological relevant MET mutants are available as recombinant human active protein kinases and for compound testing services (Figure 1; see also reverse side).



Side-by-side Comparison of MET wt and MET Mutants

All MET variants were compared side-by-side with respect to inhibition by reference inhibitor PHA665752 (Figure 2). IC₅₀ determinations of PHA665752 were performed at corresponding app. ATP K_m of each MET variant (see Table 1) in our radiometric ³³PanQinase™ Assay.

Figure 2: Differential inhibition of 10 MET variants by MET inhibitor PHA665752 at app. ATP K_m (n=2).



References

- ¹ c-Met Inhibitors with Novel Binding Mode Show Activity against Several Hereditary Papillary Renal Cell Carcinoma-related Mutations: Steven F. Bellon et al.; JBC 283, 2675-2683 (2008)
- ² The Met kinase inhibitor SU11274 exhibits a selective inhibition pattern toward different receptor mutated variants: Sylvie Berthou et al., Oncogene 23, 5387-5393 (2004)