

Kinase mutant panel performed with $^{33}\text{PanQinase}^{\text{TM}}$ assay technology.

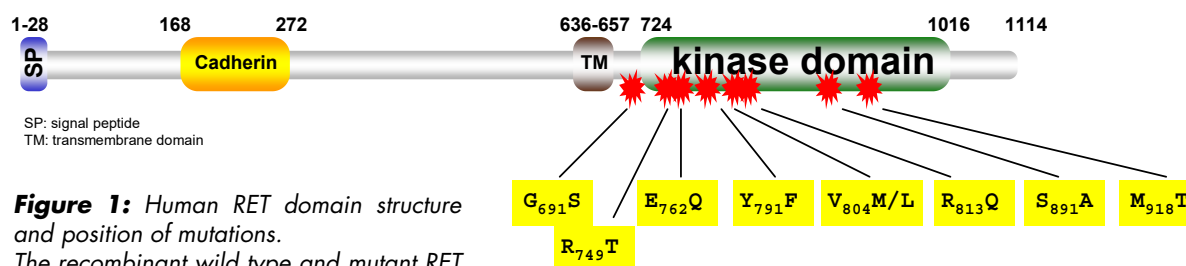
RET Kinase H658-S1114 Mutant Panel

Ret proto-oncogene

Synonyms: CDHF12, CDHR16, HSCR1, MEN2A, MEN2B, MTC1, PTC, RET51, RET-ELE1

RET is a receptor tyrosine kinase which is the catalytically active part of the signaling complex for the glial cell line-derived neurotrophic factor family of ligands. Gain-of-function mutations in RET are implicated in human cancers¹. Several physiological RET mutations have been described, RET and its mutants seem to be promising oncological targets for small molecule kinase inhibitor-based therapies².

Wild type RET as well as nine pathophysiologically relevant RET mutants are available as recombinant human active protein kinases and for compound testing services (Figure 1).

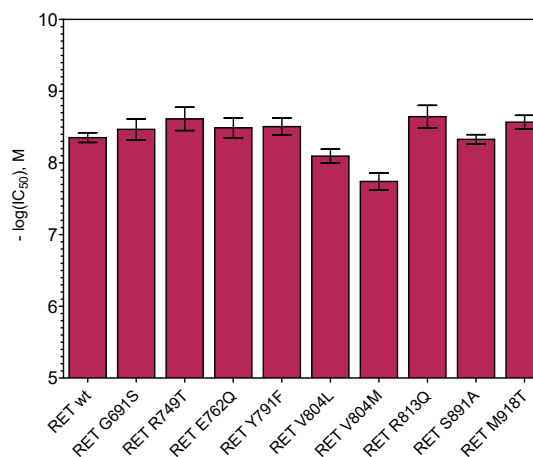


Side-by-side comparison of wild type RET and RET mutants

Wild type RET and nine different RET mutants were compared side-by-side with respect to inhibition by the reference inhibitor BIBF1120 (Vargatef) (Figure 2). IC₅₀ determinations of BIBF1120 (Vargatef) were performed at corresponding app. ATP-K_m (see Table 1) using our radiometric $^{33}\text{PanQinase}^{\text{TM}}$.

Figure 2:

Differential inhibition of 10 RET variants by the inhibitor BIBF1120 (Vargatef) at app. ATP K_m (n= 2).



References

- ¹ Structure and chemical inhibition of the RET tyrosine kinase domain: P. P. Knowles et al., J. Biol. Chem. 281, 33577-33587 (2006)
- ² Disease associated mutations at valine 804 in the RET receptor tyrosine kinase confer resistance to selective kinase inhibitors, F. Carlomagno et al., Oncogene 23, 6056-6063 (2004)