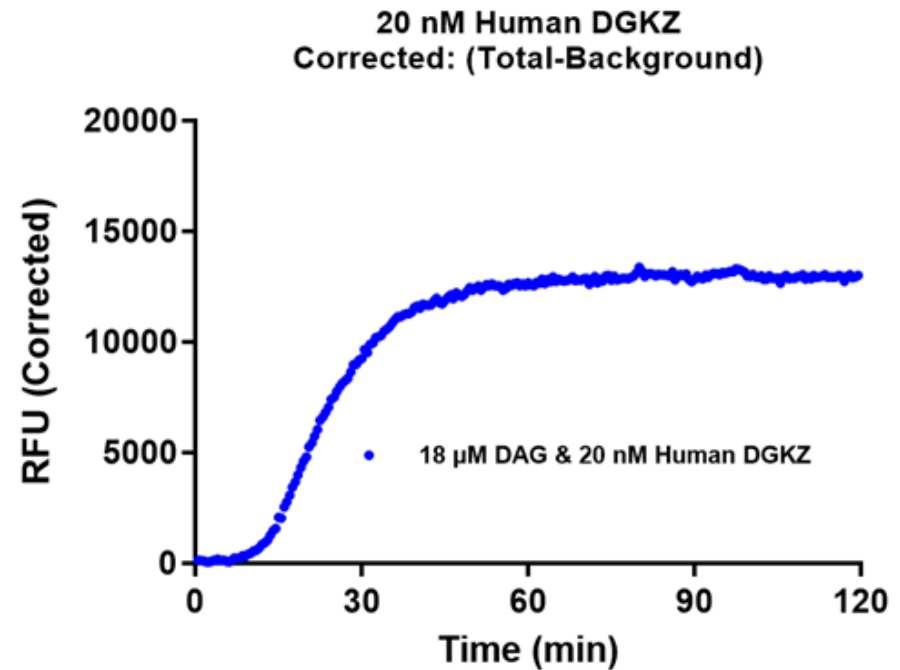
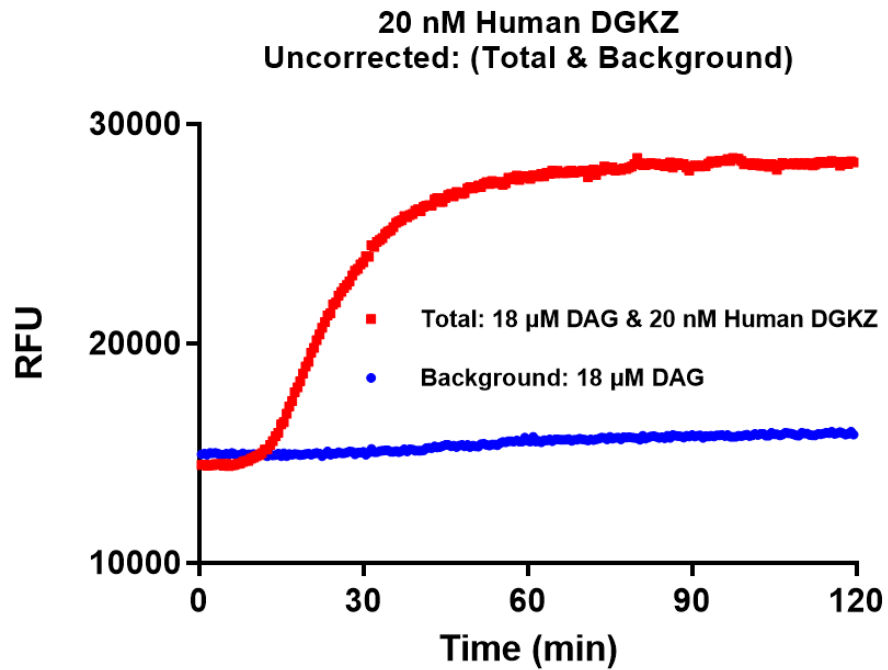


Kinase BTN-DGKζ(DGKZ)
Cat.# 12-410-20N
Substrate PhosphoSens® DAG Substrate (Assay Quant Technologies Inc.)



DAG Reaction rate: 575 ± 7 RFU/pmole/min