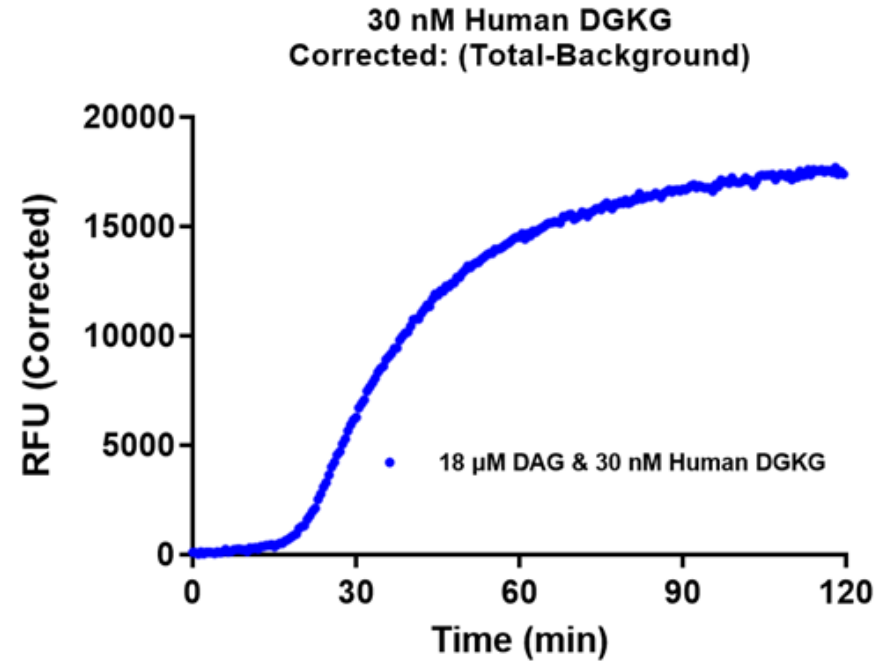
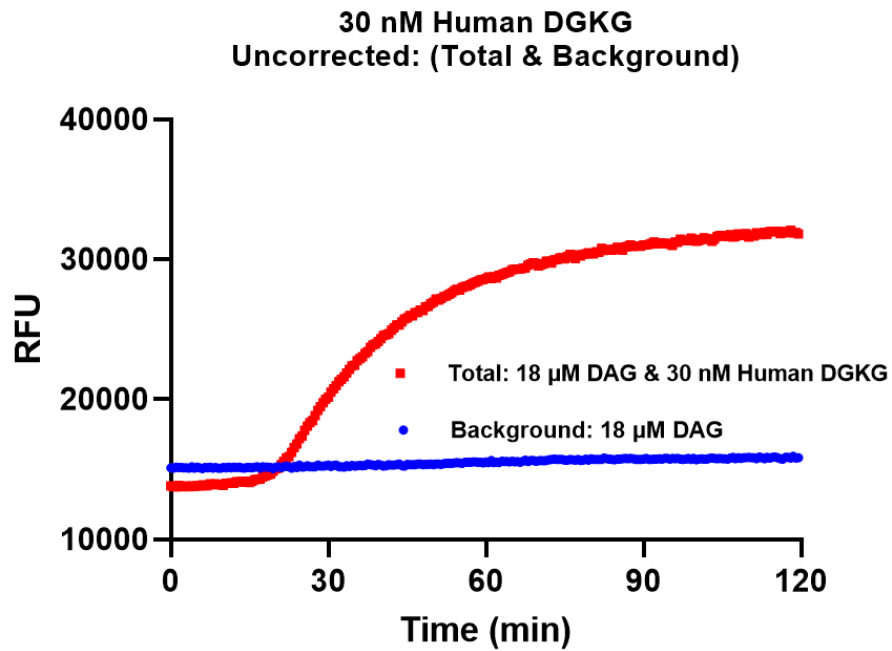


Kinase BTN-DGK γ (DGKG)
Cat.# 12-403-20N
Substrate PhosphoSens[®] DAG Substrate (Assay Quant Technologies Inc.)



DAG Reaction rate: 352 ± 4 RFU/pmole/min