

# Why should you use HTRF™?

## Are you interested in biology?

YES

NO

Are you looking for robust data?

NO

Do you want to save time?

Do you want to simplify your life in the lab?

YES

NO

YES

NO

YES

Check out [revvity.com](http://revvity.com)

Are you tired of Western Blot?

NO

Are you curious?

YES

YES

NO

Take a break


 **Go for HTRF !**

You need this!

Are you over-booked?

YES

NO

 Do you have time?

Of course you do!


NO

YES

Are you working on GPCR, PPI, cell signaling or other cutting-edge research areas?

YES

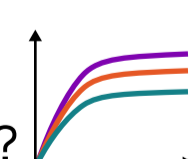
NO

 Do you want to amaze your Boss?

YES

NO

Are you interested in enzyme kinetics?



revvity