

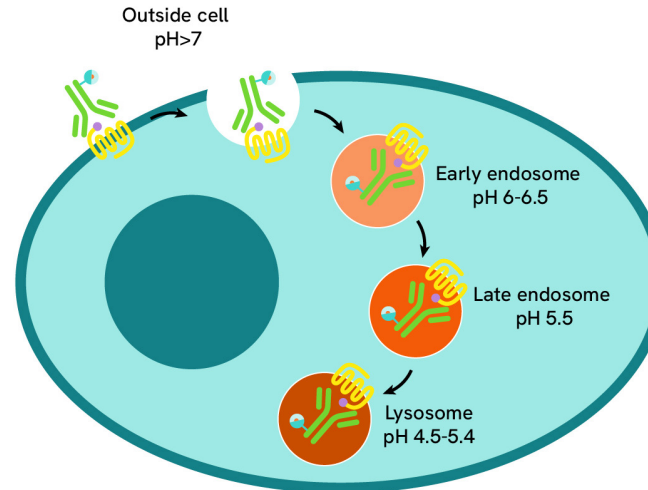
GPCR internalization: simplified detection, clear results.

Your high-throughput solution to accelerate your GPCR research is here!

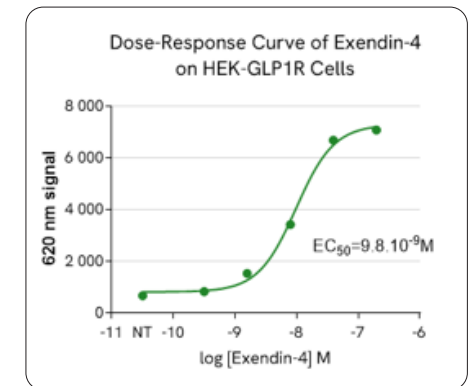
pHSense™ probes are plate reader-compatible reagents with time-resolved fluorescence (TRF) readout specifically developed for internalization studies. Fully compatible with our plate reader platforms, this technology reinforces Revvity's commitment to advancing drug discovery workflows. pHSense offers:

- Higher throughput plate-based pH sensitive internalization reagents
- Compatibility with standard plate readers
- Superior sensitivity at low endogenous expression levels

pHSense products	Part number - 96 wells
pHSense Eu Fab Anti-Human IgG	81HUEU1AA
pHSense Eu Fab Anti-Mouse IgG1	81MO1EU1AA
pHSense Eu Fab Anti-Mouse IgG2	81MO2EU1AA
pHSense Eu Anti-FLAG	81FL1EU1AA
pHSense Eu Anti-HA	81HAEU1AA
pHSense Eu SNAP Labeling Reagent	81SNEU1AA
pHSense Eu Reader Control Kit	82RCLEUPEA



While the GPCR remains at the membrane, the pHSense probe is at extracellular pH and exhibits minimal fluorescence. Upon GPCR internalization, the endosomal/lysosomal pH acidifies around the probe, which emits a strong long-lived fluorescent signal



Agonist-induced internalization of GLP1R in HEK cells