

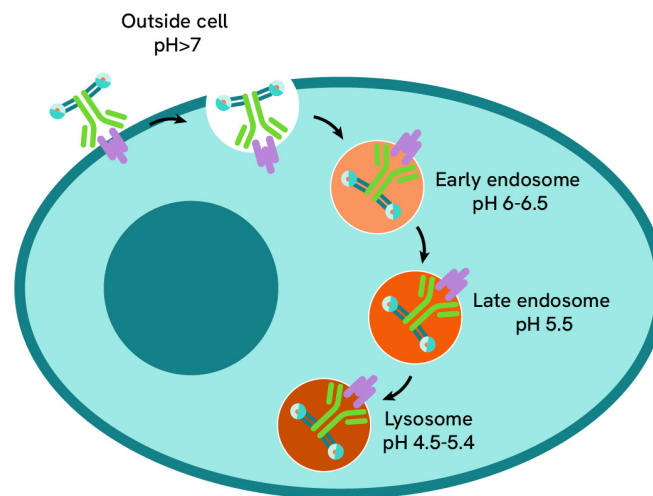
Detecting ADC internalization: from complexity to clarity.

Accelerate your ADC research with simplified detection solutions.

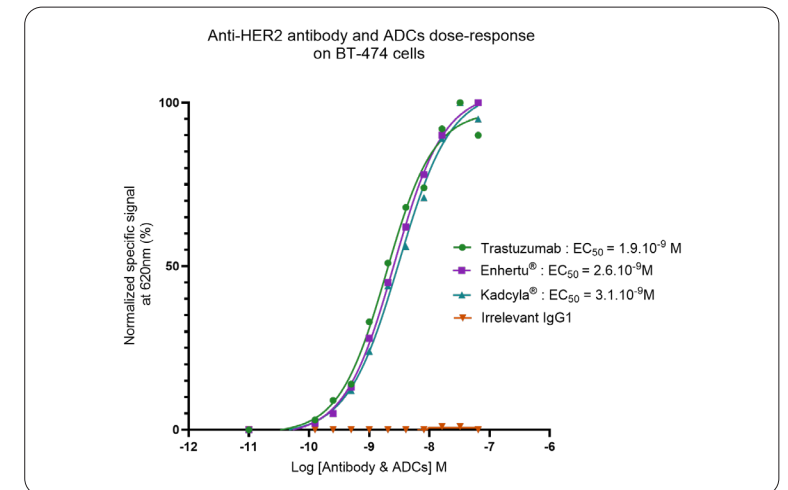
pHSense™ probes are plate reader-compatible reagents with time-resolved fluorescence (TRF) readout specifically developed for internalization studies. Fully compatible with our plate reader platforms, this technology reinforces Revvity's commitment to advancing drug discovery workflows. pHSense offers:

- Higher throughput plate-based pH sensitive internalization reagents
- Superior signal-to-background for more reliable pharmacological data
- Versatility across suspension & adherent cells

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 Reader Control Kit	82RCLEUPEA



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



Anti-HER2 internalizing antibody, Trastuzumab, and Trastuzumab-based ADCs (Kadcylla® and Enhertu®)