

## In Vivo Imaging Featured Probes Back IVISense

Product Number:      NEV20011

Storage Conditions: 2 to 8°C

Kit Components:

Component	Description	Quantity
NEV10003	IVISense Pan Cathepsin 680	1
NEV10040	IVISense Folate Receptor 680	1
NEV10054EX	IVISense Vascular 680	1
NEV10168	IVISense MMP 750 FAST	1
NEV10873	IVISense Integrin Receptor 750	1
NEV11118	IVISense 680 NHS Fluor Labeling k	1

Refer to individual components for more technical information

The information provided in this document is for reference purposes only and may not be all-inclusive. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume liability for the accuracy or completeness of the information contained herein. Users should exercise caution when handling materials as they may present unknown hazards. Revvity shall not be liable for any damages or losses resulting from handling or contact with the product, as Revvity cannot control actual methods, volumes, or conditions of use. Users are responsible for ensuring the product's suitability for their specific application. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS, OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF  
 NEV11105-R Rev01