

## Arthritis Starter Pack Option 1 IVISense

Product Number: NEV20003

Storage Conditions: 2 to 8°C

Kit Components:

Component	Description	Quantity
NEV10003	IVISense Pan Cathepsin 680	1
NEV10011EX	IVISense Vascular 750	1
NEV10020EX	IVISense Osteo 680	1
NEV10126	IVISense MMP 680	1
NEV11000	IVISense Cat K 680 FAST	1
NEV11098	IVISense Cat B 750 FAST	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