



# Great Scott!!! New brand, old prices that'll make you feel like you're back in time.

Go back in time for a limited time with IVIS® price savings.



No need for a flux capacitor; just click the button below and travel back to the future today!

Upgrade to a new and improved *in vivo* imaging instrument at introductory prices. As we complete our transformation to becoming Revvity, we're embracing our identity as a completely new company. We are thrilled to introduce our new brand to the world, and we have big plans for the future. However, we also want to celebrate our loyal customers throughout the years with this throwback savings on new IVIS® optical imaging systems. Included in this promotion are the IVIS® Lumina III Series & IVIS® Spectrum Series. Please click below for further details or reach out to your local representative.



Scan to explore *in vivo*  
imaging solutions.

Terms & Conditions: For a limited time, buy a IVIS® Lumina Series III, IVIS® Lumina LT Series III, IVIS® Lumina Series III, IVIS® Lumina XRMS, IVIS® Spectrum, or IVIS® SpectrumCT, and receive a 33% discount off the purchase price of the instrument. This promotion is open to customers in the United States. Discount will apply to purchase orders placed on or before December 31, 2023 or until promotional supplies are depleted, whichever comes first. Discount is valid toward the purchase of a new qualifying IVIS® Lumina Series III, IVIS® Lumina LT Series III, IVIS® Lumina Series III, IVIS® Lumina XRMS, IVIS® Spectrum, or IVIS® SpectrumCT. For clarity, discount is only valid on a new instrument order. Purchases of used, refurbished or demonstration instruments do not qualify for this discount. Discounts cannot be applied to existing, pending or prior orders. This promotion may not be combined with any other discounts, offers or promotions of any kind, including without limitation, other instrument price discounts, volume purchase agreements, university programs, standing order discounts or other contract pricing. Customer is responsible for any applicable charges, taxes, shipping and other related expenses; no discount will be applied to any such costs. This discount may be discontinued, suspended or amended by Revvity in its sole discretion at any time without further notice. No cash or cash equivalent. No substitutions. Offer void where prohibited, licensed, or restricted by federal, state, provincial, or local laws or regulation or agency/institutional policy. Other limitations and restrictions may apply. For details, please contact your sales representative.



| IVIS® Lumina III Series

| IVIS® Spectrum or  
Spectrum CT

Click to order