

Accessories for *in vivo* imaging.



Below is a list of a few key accessories to improve and accelerate your *in vivo* imaging studies:

Streamline *in vivo* imaging with unique accessories designed for the IVIS® imaging platform



High throughput imaging kit

Part # CLS148874

Accelerate imaging workflow with the High Throughput Imaging Kit. This kit contains two SMT-5 Smart Mouse Trays and Benchtop Posing Station, with a warming tray to keep the animals at a comfortable physiological temperature that serves as the conduit for connection of the SMT-5 tray to the RAS-4 Rodent Anesthesia System (sold separately). This kit allows for posing and queuing of up to five mice on the benchtop before placing them in the IVIS imaging system.



SRM-5 small rodent manifold

Part # CLS148872

SRM-5 small rodent manifold seamlessly delivers anesthesia to either five mice or two rats when used in conjunction with the Magnetic Quick Connect (sold separately). A robust design requires no nose cones and offers superior scavenging for reduced anesthesia gas exposure. This system provides the ultimate in flexibility and ease of use as multiple accessories can be easily removed or changed with the quick connect functionality.

Efficient rodent anesthesia system with active scavenging capabilities



RAS-4 rodent anesthesia system

Part # CLS146737

The RAS-4 Rodent Anesthesia System enables real-time *in vivo* imaging to monitor and record molecular and genetic activity in mice and rats using isoflurane gas. The RAS-4 Rodent Anesthesia System is compatible with Revvity's suite of *in vivo* imaging systems. The system facilitates efficient gas delivery via four delivery circuits while minimizing excess gas exposure to lab personnel with active scavenging capabilities.

360 Degree surface mapping



MVI-2 multi-view imager kit

Part # CLS143807

The MVI-2 Multi-view imager enables complete 360 degree 2D surface mapping of two mice simultaneously for the identification, localization, and quantification of whole body imaging of probes.

Please visit our website for a complete list of accessories <https://www.revvity.com>.

The Revvity logo, consisting of the word "revvity" in a lowercase, sans-serif font.