

**Rabbit Serum
(Normal)
Code X0902**

Recommended use	For research use only. Not for use in diagnostic procedures. Rabbit Serum (Normal) is recommended for use as control serum or reference material for rabbit serum proteins.
Reagent provided	A pool of undiluted whole sera from healthy rabbits (Danish Whites) provided in liquid form, and containing 15 mmol/L NaN ₃ . <u>Protein concentration g/L:</u> See label on vial. The protein concentration has been measured by spectrophotometry using extinction coefficient $E_{278\text{nm},1\text{cm}}^{0.1\%} = 1.00$ A fine precipitate of serum proteins may form in the serum. This precipitate may be removed, if required, by passing the serum through a 0.2 µm filter immediately before use.
Precautions	<ol style="list-style-type: none"> 1. The device is not intended for clinical use including diagnosis, prognosis, and monitoring of a disease state, and it must not be used in conjunction with patient records or treatment. 2. This product contains sodium azide (NaN₃), a chemical highly toxic in pure form. At product concentrations, though not classified as hazardous, sodium azide may react with lead and copper plumbing to form highly explosive build-ups of metal azides. Upon disposal, flush with large volumes of water to prevent metal azide build-up in plumbing. 3. As with any product derived from biological sources, proper handling procedures should be used. 4. Wear appropriate Personal Protective Equipment to avoid contact with eyes and skin. 5. Unused solution should be disposed of according to local, State and Federal regulations.
Storage	Store at 2-8 °C. Do not use after expiration date stamped on vial. If unexpected staining is observed which cannot be explained by variations in laboratory procedures and a problem with the product is suspected, contact our Technical Services.

Explanation of symbols

 REF	Catalogue number	 LOT	Batch code
	Consult instructions for use		Use by
 2°C-8°C	Temperature limitation		Manufacturer