

Mouse IgG2b Isotype Control (ATTO 488 conjugate)

This antibody is covered by our [Worry-Free Guarantee](#).

Ordering Information

[Order Online »](#)

ALX-804-877TD-C050	ATTO 488 - 50µg
--------------------	-----------------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored as recommended.
Handling	Do not freeze. Protect from light.
Short Term Storage	+4°C
Long Term Storage	+4°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application Notes	Negative control in assays where mouse IgG2b (ATTO 488) antibodies are used.
Clone	A-1
Formulation	Liquid. In PBS containing 0.02% sodium azide.
Host	Mouse
Isotype	IgG2b κ
Species Reactivity	Species independent
Technical Info / Product Notes	ATTO-fluorescent antibodies show increased photostability, outstanding brightness and intense signals. ATTO dyes are thermally stable, resistant to environmental changes and show no significant isomerization. ATTO 488 shows bright green fluorescence (λ_{abs} (max): 501nm, λ_{em} (max): 523nm, ϵ_{max} : 90'000).

Worry-free Guarantee

This antibody is covered by our [Worry-Free Guarantee](#)