



EAGLE
BIOSCIENCES

Monoclonal Mouse Anti-Human BRAF V600E Antibody

Catalog Number:

IHC600-100 0.1 mL, Concentrate

IHC600-1 1 mL, Concentrate

IHC600-7 7 mL, Predilute

For Research Use Only. Not for use in diagnostic procedures.

v. 1.0 08/12/2021

Eagle Biosciences, Inc.

20A Northwest Blvd., Suite 112, Nashua, NH 03063

Phone: 617-419-2019 Fax: 617-419-1110

www.EagleBio.com

