

**MicroTools™ Starter Kit 2**

MicroTools Kit 2 on 25 mm rods, preloaded into reusable B1A magnetic bases

Cat. No.	Amount
T2SK-1	1 Kit

**For general laboratory use.****Shipping:** shipped at ambient temperature**Storage Conditions:** store at ambient temperature**Shelf Life:** n/a**Content:**

- 20 Individual MicroTools on 25 mm rods, preloaded into B1A-R magnetic bases
- Magnetic Push-button Wand (for easily switching between tools)
- Serrated-end tweezers
- Holder and case
- Tool location placard

**Related Products:**MicroTools™ Starter Kit 1, #T1SK-1  
MicroTools™ Starter Kit 3, #T3SK-1**nonsaleable to USA and Japan**