

**MicroTools™ Starter Kit 1**

MicroTools Kit 1 on 25 mm rods, preloaded into reusable B1A magnetic bases

Cat. No.	Amount
T1SK-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 2, #T2SK-1

MicroTools™ Starter Kit 3, #T3SK-1

nonsaleable to USA and Japan