



## 2'-Fluoro-dNTP Bundle

(2'F-dNTP Bundle)

2'-F-2'-dNTP

2'-Fluoro-2'-deoxynucleoside-5'-triphosphates, Sodium salts

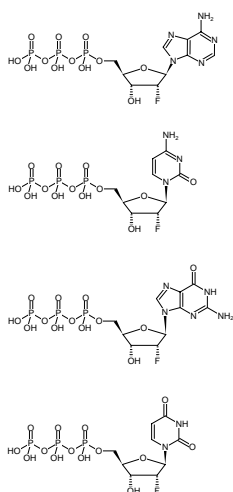
2'-Deoxy-2'-fluoronucleoside-5'-triphosphates

Cat. No.	Amount
NU-10405-20	4 x 50 µl (4 x 5 µmol)

### Description:

2'-Fluoro-dNTP Bundle is a set of 2'-Fluoro-2'-deoxynucleoside-5'-triphosphates with 5 µmol (50 µl/100mM) of each.

2'Fluoro-dATP (#NU-151)  
 2'Fluoro-dCTP (#NU-1214)  
 2'Fluoro-dGTP (#NU-1216)  
 2'Fluoro-dUTP (#NU-1215)



Structural formula of 2'-Fluoro-dNTP Bundle

**For *in vitro* use only!**

**Shipping:** shipped on blue ice

**Storage Conditions:** store at -20 °C

**Additional Storage Conditions:** store dark and dry

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

**Shelf Life:** 12 months after date of delivery

### Molecular Formula:

2'Fluoro-dATP: C<sub>10</sub>H<sub>15</sub>FN<sub>5</sub>O<sub>12</sub>P<sub>3</sub> (free acid)

2'Fluoro-dCTP: C<sub>9</sub>H<sub>15</sub>FN<sub>3</sub>O<sub>13</sub>P<sub>3</sub> (free acid)

2'Fluoro-dGTP: C<sub>10</sub>H<sub>15</sub>FN<sub>5</sub>O<sub>13</sub>P<sub>3</sub> (free acid)

2'Fluoro-dUTP: C<sub>9</sub>H<sub>14</sub>FN<sub>2</sub>O<sub>14</sub>P<sub>3</sub> (free acid)

### Molecular Weight:

2'Fluoro-dATP: 509.17 g/mol (free acid)

2'Fluoro-dCTP: 485.15 g/mol (free acid)

2'Fluoro-dGTP: 525.17 g/mol (free acid)

2'Fluoro-dUTP: 486.13 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

**Form:** clear aqueous solution

**Concentration:** 100 mM