

## Incucyte® 实时活细胞分析系统 试剂与耗材

活细胞分析系统搭配 Incucyte® 试剂和耗材，为广泛的细胞健康、运动、功能等应用提供一体化解决方案。

Incucyte® 试剂不仅能够提高效率、保持细胞健康，还不会对细胞的增殖和形态产生干扰。致力于提供满足您应用需求的解决方案！

### 细胞凋亡与细胞毒性

产品	颜色	数量	兼容性	目录号
Incucyte® Annexin V 绿色染料		1 支: 100 tests	SX1, S3, SX5, CX3	4642
Incucyte® Annexin V 红色染料		1 支: 100 tests	SX1, S3, SX5 (配置绿色 / 红色光学模块)	4641
Incucyte® Annexin V NIR 染料		1 支: 100 tests	SX5, CX3	4768
Incucyte® Annexin V 橙色染料		1 支: 100 tests	SX5, CX3	4759
Incucyte® Caspase-3/7 绿色染料		1 支: 20 µL	SX1, S3, SX5, CX3	4440
Incucyte® Caspase-3/7 红色染料		1 支: 20 µL	SX1, S3, SX5 (配置绿色 / 红色光学模块)	4704
Incucyte® Cytotox 绿色染料		5 支: 每支 5 µL	SX1, S3, SX5, CX3	4633
Incucyte® Cytotox 红色染料		5 支: 每支 5 µL	SX1, S3, SX5 (配置绿色 / 红色光学模块)	4632
Incucyte® Cytotox NIR 染料		1 支: 100 µL	SX5, CX3	4846

### 细胞增殖

产品	颜色	数量	兼容性	目录号
Incucyte® Nuclight 红色慢病毒 2.0 试剂 (puro)		一支: 0.6 mL	SX1, S3, SX5 (配置绿色 / 红色光学模块)	BA-04887
Incucyte® Nuclight 绿色慢病毒 2.0 试剂 (puro)		一支: 0.6 mL	SX1, S3, SX5, CX3	BA-04888
Incucyte® Nuclight 红色慢病毒 2.0 试剂 (bleo)		一支: 0.6 mL	SX1, S3, SX5 (配置绿色 / 红色光学模块)	BA-04889
Incucyte® Nuclight 绿色慢病毒 2.0 试剂 (bleo)		一支: 0.6 mL	SX1, S3, SX5, CX3	BA-04890
Incucyte® Nuclight 橙色慢病毒 2.0 试剂 (puro)		一支: 0.6 mL	SX5, CX3	BA-04891
Incucyte® NIR 慢病毒试剂 (puro)		一支: 0.2 mL	SX5, CX3	4805
Incucyte® Nuclight 快速红色染料		1 支: 50 µL	SX1, S3, SX5 (配置绿色 / 红色光学模块)	4717
Incucyte® Nuclight 快速 NIR 染料		1 支: 50 µL	SX5, CX3	4804

### 细胞周期

产品	颜色	数量	兼容性	目录号
Incucyte® 细胞周期绿 / 红慢病毒 2.0 试剂		1 支: 0.2 mL	SX1, S3, SX5 (配置绿色 / 红色光学模块)	BA-04892
Incucyte® 细胞周期绿 / 橙慢病毒 2.0 试剂		1 支: 0.2 mL	SX5, CX3	BA-04889

## 免疫细胞杀伤、激活和增殖

产品	颜色	数量	兼容性	目录号
Incucyte® Annexin V 绿色染料		1 支: 100 tests	SX1, S3, SX5, CX3	4642
Incucyte® Annexin V 红色染料		1 支: 100 tests	SX1, S3, SX5 (配置绿色 / 红色光学模块)	4641
Incucyte® Annexin V NIR 染料		1 支: 100 tests	SX5, CX3	4768
Incucyte® Annexin V 橙色染料		1 支: 100 tests	SX5, CX3	4759
Incucyte® Caspase-3/7 绿色染料		1 支: 20 µL	SX1, S3, SX5, CX3	4440
Incucyte® Caspase-3/7 红色染料		1 支: 20 µL	SX1, S3, SX5 (配置绿色 / 红色光学模块)	4704
Incucyte® Cytolight 快速绿色染料		1 支: 15 µg	SX1, S3, SX5, CX3	4705
Incucyte® Cytolight 快速红色染料		5 支: 50 µg	SX1, S3, SX5 (配置绿色 / 红色光学模块)	4706
Incucyte® Cytolight 快速橙色染料		1 支: 1 mg	SX5, CX3	4839

## 划痕细胞迁移和侵袭

产品	数量	兼容性	目录号
Incucyte® Imagelock 96 孔板	50 块板	SX1, S3, SX5, CX3	BA-04857
	10 块板		BA-04856
Incucyte® 96 孔划痕器	1 个	SX1, S3, SX5, CX3	4563
Incucyte® 细胞迁移组合包	1 个	SX1, S3, SX5, CX3	BA-04858

## 趋化性细胞迁移和侵袭

产品	数量	兼容性	目录号
Incucyte® Clearview 96 孔趋化板	1 块板	SX1, S3, SX5, CX3	4582
	10 块板		4648
	50 块板		4599
Incucyte® ClearView 96 孔储板	10 块板	SX1, S3, SX5, CX3	4600
	100 块板		4601

## 免疫细胞化学

产品	颜色	数量	兼容性	目录号
Incucyte® 小鼠 IgG1 Fabfluor-488 抗体标记试剂盒 ■ Incucyte® 小鼠 IgG1 FabFluor-488 染料 ■ Incucyte® Opti-Green 背景抑制剂		1 个	SX1, S3, SX5, CX3	4745
Incucyte® 小鼠 IgG2a Fabfluor-488 抗体标记试剂盒 ■ Incucyte® 小鼠 IgG12a FabFluor-488 染料 ■ Incucyte® Opti-Green 背景抑制剂		1 个	SX1, S3, SX5, CX3	4743
Incucyte® 小鼠 IgG2b Fabfluor-488 抗体标记试剂盒 ■ Incucyte® 小鼠 IgG12b FabFluor-488 染料 ■ Incucyte® Opti-Green 背景抑制剂		1 个	SX1, S3, SX5, CX3	4744
Incucyte® 小鼠 IgG1 Fabfluor-555 抗体标记试剂盒 ■ Incucyte® 小鼠 IgG1 FabFluor-555 染料 ■ Incucyte® Opti-Orange 背景抑制剂		1 个	SX5, CX3	BA-04873
Incucyte® 小鼠 IgG1 Fabfluor-594 抗体标记染料		1 支: 50 µg	SX1, S3, SX5 (配置绿色 / 红色光学模块)	4844
Incucyte® 小鼠 IgG2a Fabfluor-594 抗体标记染料		1 支: 50 µg	SX1, S3, SX5 (配置绿色 / 红色光学模块)	BA-04863

## 神经细胞活性

产品	数量	兼容性	目录号
Incucyte® Neuroburst 橙色慢病毒试剂	一支: 2 mL	SX5, CX3	4736
Incucyte® Neuroburst 橙色试剂盒	1 个	SX5, CX3	4761

## 抗体内化

产品	颜色	数量	兼容性	目录号
Incucyte® 人 Fabfluor-pH 红色抗体标记染料	■	一支: 50 µg	SX1, S3, SX5 (配置绿色 / 红色光学模块)	4722
Incucyte® 人 Fabfluor-pH 橙色抗体标记染料	■	一支: 50 µg	SX5, CX3	4812
Incucyte® 小鼠 IgG1Fabfluor-pH 红色抗体标记染料	■	一支: 50 µg	SX1, S3, SX5 (配置绿色 / 红色光学模块)	4723
Incucyte® 小鼠 IgG1Fabfluor-pH 橙色抗体标记染料	■	一支: 50 µg	SX5, CX3	BA-04885
Incucyte® 小鼠 IgG2a Fabfluor-pH 红色抗体标记染料	■	一支: 50 µg	SX1, S3, SX5 (配置绿色 / 红色光学模块)	4750
Incucyte® 小鼠 IgG2a Fabfluor-pH 橙色抗体标记染料	■	一支: 50 µg	SX5, CX3	BA-04886
Incucyte® 小鼠 IgG2b Fabfluor-pH 红色抗体标记染料	■	一支: 50 µg	SX1, S3, SX5 (配置绿色 / 红色光学模块)	4751
Incucyte® 大鼠 Fabfluor-pH 红色抗体标记染料	■	一支: 50 µg	SX1, S3, SX5 (配置绿色 / 红色光学模块)	4737

## 细胞代谢

产品	颜色	数量	兼容性	目录号
Incucyte® CytoATP 慢病毒试剂盒 ■ Incucyte® CytoATP 慢病毒标记试剂 ■ Incucyte® CytoATP Non-binding Control (puro)	■	1 个	SX5 (配置 SX5 代谢光学模块)	4772

## 吞噬作用

产品	颜色	数量	兼容性	目录号
pHrodo® 红色细胞标记试剂盒	■	1 个	SX1, S3, SX5 (配置绿色 / 红色光学模块)	4649
pHrodo® 橙色细胞标记试剂盒	■	1 个	SX5, CX3	4766
Incucyte® pHrodo® 大肠杆菌红色荧光颗粒	■	一支: 2 mg	SX1, S3, SX5 (配置绿色 / 红色光学模块)	4615
Incucyte® pHrodo® 大肠杆菌绿色荧光颗粒	■	一支: 2 mg	SX1, S3, SX5, CX3	4616

## 细胞外囊泡 (EV)

产品	数量	兼容性	目录号
Incucyte® 细胞外囊泡 (EV) 绿色标记试剂盒	1 个	SX1, S3, SX5, CX3	BA-04877

# 联系我们

更多联系信息，请访问

[www.sartorius.com.cn](http://www.sartorius.com.cn)

赛多利斯莱珀思（上海）贸易有限公司

邮箱 [leadscn@sartorius.com](mailto:leadscn@sartorius.com)

服务热线 400 920 9889 | 800 820 9889

