

推荐货号	规格	产品名称	代谢分类	所享折扣
ab83428	100 test	Oxaloacetate Assay Kit	中间代谢	五折
ab102528	100 test	Isocitrate Dehydrogenase Assay Kit (Colorimetric)	中间代谢	五折
ab183305	100 test	Malate Dehydrogenase Activity Assay Kit (Colorimetric)	中间代谢	五折
ab219926	200 test	Glutathione Peroxidase Activity Assay Kit (Fluorometric)	抗氧化	五折
ab197242	96 test	Extracellular Oxygen Consumption Reagent	氧化应激	五折
ab197243	4 x 96 tests	Extracellular Oxygen Consumption Assay	氧化应激	五折
ab232857	1 x 96 tests	Mitochondrial Stress Test Complete Assay Kit	氧化应激	五折
ab65347	100 test	L-Amino Acid Assay Kit	氨基酸代谢	五折
ab83394	100 test	L-Alanine Assay Kit	氨基酸代谢	五折
ab102512	100 test	Aspartate Assay Kit	氨基酸代谢	五折
ab126286	100 test	Inosine Quantification Assay Kit (Fluorometric)	氨基酸代谢	五折
ab204695	100 test	Adenosine Deaminase (ADA) Activity Assay Kit (Fluorometric)	氨基酸代谢	五折
ab204707	100 test	Purine Nucleoside Phosphorylase Activity Assay Kit (Colorimetric)	氨基酸代谢	五折
ab102532	100 test	Glucose Dehydrogenase Assay Kit (Colorimetric)	糖代谢	五折
ab133118	96 test	Sphingomyelin Assay Kit	细胞信号通路相关	五折
ab183306	100 test	Phospholipase D Activity Assay Kit (Colorimetric)	细胞信号通路相关	五折
ab190554	200 test	Acidic Sphingomyelinase Assay Kit (Fluorometric)	细胞信号通路相关	五折
ab103967	300 µl	Triglyceride Standard	脂代谢	五折
ab105138	10 ml	2X LDL/VLDL Precipitation Buffer	脂代谢	五折
ab83355	100 test	ATP Assay Kit (Colorimetric/Fluorometric)	中间代谢	六折
ab83359	100 test	ADP Assay Kit (Colorimetric/Fluorometric)	中间代谢	六折
ab83391	100 test	Malate Assay Kit	中间代谢	六折
ab83431	100 test	Alpha Ketoglutarate (alpha KG) Assay Kit	中间代谢	六折
ab83432	100 test	Pyruvate Kinase Assay Kit	中间代谢	六折
ab102516	100 test	Fumarate Assay Kit	中间代谢	六折
ab102527	100 test	Glutamate Dehydrogenase Activity Assay Kit	中间代谢	六折
ab102533	100 test	Alcohol Dehydrogenase Assay Kit	中间代谢	六折
ab113849	300 test	Luminescent ATP Detection Assay Kit	中间代谢	六折
ab174094	100 test	Glycerol-3-phosphate (G3P) Assay Kit (Colorimetric)	中间代谢	六折
ab196994	100 test	Aldolase Activity Assay Kit (Colorimetric)	中间代谢	六折
ab197001	100 test	Triose Phosphate Isomerase (TPI) Activity Assay Kit (Colorimetric)	中间代谢	六折
ab204718	100 test	Succinate Assay Kit (Colorimetric)	中间代谢	六折
ab83396	100 test	Citrate Assay Kit	中间代谢	六折
ab110671	1 x 96 tests	Pyruvate dehydrogenase (PDH) Combo (Activity + Profiling) Microplate Assay Kit	中间代谢	六折
ab140361	1 x 96 tests	Lactate Dehydrogenase B / LDH-B Activity Assay	中间代谢	六折
ab176722	200 test	Glucose 6 Phosphate Dehydrogenase Activity Assay Kit (Fluorometric)	中间代谢	六折
ab196992	100 test	Fumarase Activity Assay Kit (Colorimetric)	中间代谢	六折
ab211071	100 test	6 Phosphogluconate Assay Kit (Colorimetric)	中间代谢	六折
ab211087	100 test	ADP Assay Kit (Fluorometric)	中间代谢	六折
ab241016	100 test	6-Phosphogluconate Dehydrogenase Assay Kit	中间代谢	六折
ab241031	100 test	Monoamine Oxidase (MAO) Assay Kit	中间代谢	六折
ab65322	100 test	Glutathione Assay Kit (Fluorometric)	抗氧化	六折

ab102530	100 test	Glutathione Peroxidase Assay Kit (Colorimetric)	抗氧化	六折
ab138881	100 test	GSH/GSSG Ratio Detection Assay Kit (Fluorometric - Green)	抗氧化	六折
ab205811	100 test	GSH/GSSG Ratio Detection Assay Kit II (Fluorometric - Green)	抗氧化	六折
ab65349	100 test	NADP/NADPH Assay Kit	氧化应激	六折
ab112132	100 test	Intracellular glutathione (GSH) Detection Assay Kit	氧化应激	六折
ab112139	500 test	Lysyl Oxidase Activity Assay Kit (Fluorometric)	氧化应激	六折
ab176724	250 test	NADP/NADPH Assay Kit (Fluorometric)	氧化应激	六折
ab186033	400 test	Total NADP and NADPH Assay Kit (Colorimetric)	氧化应激	六折
ab197244	96 test	Glycolysis Assay [Extracellular acidification]	氧化应激	六折
ab197245	96 test	Intracellular Oxygen Concentration Assay	氧化应激	六折
ab65339	100 test	Creatine Assay Kit (Colorimetric/Fluorometric)	氨基酸代谢	六折
ab65340	100 test	Creatinine Assay Kit	氨基酸代谢	六折
ab65344	100 test	Uric Acid Assay Kit	氨基酸代谢	六折
ab83360	100 test	Ammonia Assay Kit	氨基酸代谢	六折
ab83362	100 test	Urea Assay Kit	氨基酸代谢	六折
ab83374	100 test	Branched Chain Amino Acid Assay Kit / BCAA Assay Kit	氨基酸代谢	六折
ab83376	100 test	Phenylalanine Assay Kit	氨基酸代谢	六折
ab83389	100 test	Glutamate Assay Kit	氨基酸代谢	六折
ab105134	100 test	Alanine Transaminase Activity Assay Kit (Colorimetric/Fluorometric)	氨基酸代谢	六折
ab105135	100 test	Aspartate Aminotransferase Activity Assay Kit	氨基酸代谢	六折
ab107922	100 test	Asparaginase Activity Assay Kit	氨基酸代谢	六折
ab133081	96 test	Dipeptidyl peptidase IV (DPP4) Inhibitor Screening Assay Kit	氨基酸代谢	六折
ab138883	200 test	Glutamate Assay Kit (Fluorometric)	氨基酸代谢	六折
ab138888	200 test	Neuraminidase Assay Kit (Fluorometric - Blue)	氨基酸代谢	六折
ab155901	100 test	Creatine Kinase Activity Assay Kit (Colorimetric)	氨基酸代谢	六折
ab180877	100 test	Arginase Activity Assay Kit (Colorimetric)	氨基酸代谢	六折
ab197011	100 test	Glutamine Assay Kit (Colorimetric)	氨基酸代谢	六折
ab204697	100 test	Urease Activity Assay Kit (Colorimetric)	氨基酸代谢	六折
ab211093	100 test	Adenosine Deaminase (ADA) Activity Assay Kit (Colorimetric)	氨基酸代谢	六折
ab102509	100 test	Ammonia Assay Kit - Modified Berthelot - (Colorimetric)	氨基酸代谢	六折
ab138878	1 kit	Aspartate Aminotransferase Activity Assay Kit (Fluorometric)	氨基酸代谢	六折
ab197002	100 test	Adenosylhomocysteinase (AHCY) Activity Assay Kit (Fluorometric)	氨基酸代谢	六折
ab204694	100 test	Adenosylhomocysteinase (AHCY) Inhibitor Assay Kit (Fluorometric)	氨基酸代谢	六折
ab211098	100 test	Tryptophan Assay Kit (Fluorometric)	氨基酸代谢	六折
ab211099	100 test	Cysteine Assay Kit (Fluorometric)	氨基酸代谢	六折
ab211100	100 test	Glycine Assay Kit (Fluorometric)	氨基酸代谢	六折
ab241000	100 test	Phenylalanine Assay Kit (Colorimetric)	氨基酸代谢	六折
ab252893	100 test	Glutamate Assay Kit (Fluorometric)	氨基酸代谢	六折
ab65330	100 test	L-Lactate Assay Kit (Colorimetric/Fluorometric)	糖代谢	六折
ab65331	100 test	L-Lactate Assay Kit (Colorimetric)	糖代谢	六折
ab65333	100 test	Glucose Assay Kit	糖代谢	六折
ab65342	100 test	Pyruvate Assay Kit	糖代谢	六折
ab65343	100 test	Ethanol Assay Kit	糖代谢	六折

ab65620	100 test	Glycogen Assay Kit	糖代谢	六折
ab83375	100 test	Sialic Acid (NANA) Assay Kit	糖代谢	六折
ab83380	100 test	Fructose Assay Kit	糖代谢	六折
ab83382	100 test	Galactose Assay Kit	糖代谢	六折
ab83429	100 test	D-Lactate Assay Kit (Colorimetric)	糖代谢	六折
ab102517	100 test	Glucose Assay Kit - reducing agent compatible	糖代谢	六折
ab102526	500 test	LDH Assay Kit / Lactate Dehydrogenase Assay Kit (Colorimetric)	糖代谢	六折
ab102529	100 test	Glucose 6 Phosphate Dehydrogenase Assay Kit (Colorimetric)	糖代谢	六折
ab138884	500 test	Glucose Oxidase Assay Kit (Fluorometric)	糖代谢	六折
ab155898	100 test	6-Phosphofructokinase Activity Assay Kit (Colorimetric)	糖代谢	六折
ab169559	100 test	Glucose Assay Kit (Fluorometric, High Sensitivity)	糖代谢	六折
ab174093	100 test	Alpha-Glucosidase Activity Assay Kit (Colorimetric)	糖代谢	六折
ab197003	100 test	Dihydroxyacetone Phosphate (DHAP) Assay Kit (Fluorometric)	糖代谢	六折
ab204702	50 test	Glucose Uptake Assay Kit (Cell-based)	糖代谢	六折
ab204705	100 test	N-Acetylglucosaminidase (beta-NAG) Activity Assay Kit (Colorimetric)	糖代谢	六折
ab204722	100 test	Dipeptidyl peptidase IV (DPP4) Activity Assay Kit (Fluorometric)	糖代谢	六折
ab204732	100 test	Glyceraldehyde 3 Phosphate Dehydrogenase Activity Assay Kit (Colorimetric)	糖代谢	六折
ab211114	100 test	Hexokinase II Inhibitor Assay Kit (Colorimetric)	糖代谢	六折
ab235976	1 x 96 tests	2-NBDG Glucose Uptake Assay Kit	糖代谢	六折
ab239711	100 test	Cytochrome C Oxidase Assay Kit	糖代谢	六折
ab83384	100 test	Lactose Assay Kit	糖代谢	六折
ab83426	100 test	Glucose-6-Phosphate Assay Kit (Colorimetric)	糖代谢	六折
ab107923	100 test	Glucose-6-Phosphate Assay Kit - High Sensitivity (Fluorometric)	糖代谢	六折
ab118968	100 test	D-Sorbitol Assay Kit (Colorimetric)	糖代谢	六折
ab155891	100 test	Total Carbohydrate Assay Kit - Quantification	糖代谢	六折
ab155892	100 test	Glucose-1-Phosphate Assay Kit (Colorimetric)	糖代谢	六折
ab155893	100 test	ALDH Activity Assay Kit (Colorimetric)	糖代谢	六折
ab155894	100 test	ALDH Activity Assay Kit (Fluorometric)	糖代谢	六折
ab155896	100 test	Phosphoglucomutase Activity Assay Kit (Colorimetric)	糖代谢	六折
ab155897	100 test	Phosphoglucose Isomerase Activity Assay Kit (Colorimetric)	糖代谢	六折
ab169557	100 test	L-Lactate Assay Kit (Fluorometric, High Sensitivity)	糖代谢	六折
ab197000	500 test	Lactate Dehydrogenase (LDH) Assay Kit (Fluorometric)	糖代谢	六折
ab239712	100 test	Citrate Synthase Assay Kit	糖代谢	六折
ab241006	100 test	Methylglyoxal Assay Kit	糖代谢	六折
ab65345	100 test	Choline/Acetylcholine Assay Kit	细胞信号通路相关	六折
ab83377	100 test	Phosphatidylcholine Assay Kit (Colorimetric/Fluorometric)	细胞信号通路相关	六折
ab138876	200 test	Sphingomyelinase Assay Kit (Colorimetric)	细胞信号通路相关	六折
ab138877	100 test	Sphingomyelin Assay Kit (Fluorometric)	细胞信号通路相关	六折
ab241038	100 test	Lipoxygenase Assay Kit	细胞生化检测	六折
ab102523	100 test	Amylase Assay Kit (Colorimetric)	细胞生物活动相关	六折

ab133089	96 test	Secretory Phospholipase A2 Assay Kit	细胞生物活动相关	六折
ab133090	96 test	Cytosolic Phospholipase A2 Assay Kit	细胞生物活动相关	六折
ab168540	100 test	MitoBiogenesis™ Flow Cytometry Kit	细胞生物活动相关	六折
ab65336	100 test	Triglyceride Assay Kit - Quantification	脂代谢	六折
ab65337	100 test	Free Glycerol Assay Kit	脂代谢	六折
ab65341	100 test	Free Fatty Acid Assay Kit - Quantification	脂代谢	六折
ab65390	100 test	Cholesterol Assay Kit - HDL and LDL/VLDL	脂代谢	六折
ab83390	100 test	beta Hydroxybutyrate (beta HB) Assay Kit (Colorimetric)	脂代谢	六折
ab83392	100 test	L-Carnitine Assay Kit	脂代谢	六折
ab87546	100 test	Acetyl CoA Assay Kit	脂代谢	六折
ab102513	100 test	Adipogenesis Detection Assay Kit (Colorimetric/Fluorometric)	脂代谢	六折
ab102515	100 test	Cholesterol/Cholesteryl Ester Assay Kit - Detection	脂代谢	六折
ab118183	96 test	Fatty Acid Oxidation Assay Kit (flow cytometry)	脂代谢	六折
ab133102	1 kit	Adipogenesis Assay Kit (Cell-Based)	脂代谢	六折
ab133115	1 kit	Adipolysis Assay Kit	脂代谢	六折
ab133127	1 kit	LDL Uptake Assay Kit (Cell-Based)	脂代谢	六折
ab174092	100 test	Free Glycerol Assay Kit (Fluorometric, High Sensitivity)	脂代谢	六折
ab174095	100 test	Glycerol-3-Phosphate Dehydrogenase (G3PDH) Assay Kit (Colorimetric)	脂代谢	六折
ab176768	100 test	Free Fatty Acid Uptake Assay Kit (Fluorometric)	脂代谢	六折
ab178780	100 test	Triglyceride Assay Kit (Fluorometric, Reducing Samples)	脂代谢	六折
ab185433	100 test	Lipolysis Assay Kit (Colorimetric)	脂代谢	六折
ab185434	100 test	Lipolysis Assay Kit (Fluorometric)	脂代谢	六折
ab196985	100 test	Cholesterol Efflux Assay Kit (Cell-based)	脂代谢	六折
ab204701	100 test	HMG-CoA Reductase Activity Assay Kit (Colorimetric)	脂代谢	六折
ab204721	100 test	Lipoprotein Lipase Assay Kit (Fluorometric)	脂代谢	六折
ab217602	55 test	Fatty Acid Oxidation Assay	脂代谢	六折
ab222944	55 test	Fatty Acid Oxidation Complete Assay Kit	脂代谢	六折
ab234050	100 test	Phospholipid Assay Kit (Colorimetric/Fluorometric)	脂代谢	六折
ab236212	1 kit	Cholesterol Uptake Assay Kit	脂代谢	六折
ab239702	100 test	Bile Acid Assay Kit (Colorimetric)	脂代谢	六折
ab102524	100 test	Lipase Assay Kit (Colorimetric)	脂代谢	六折
ab102525	100 test	Lipase Assay Kit II (Colorimetric)	脂代谢	六折
ab118969	100 test	Lipase Assay Kit III (Fluorometric)	脂代谢	六折
ab133131	1 kit	Hepatic Lipid Accumulation/ Steatosis Assay Kit	脂代谢	六折
ab204717	100 test	HDL Uptake Assay Kit (Fluorometric)	脂代谢	六折
ab241005	100 test	Phosphatidylethanolamine Assay Kit	脂代谢	六折
ab118970	100 test	Lipid Peroxidation (MDA) Assay Kit (Colorimetric/Fluorometric)	脂肪氧化	六折
ab133085	100 test	Lipid Hydroperoxide (LPO) Assay Kit	脂肪氧化	六折
ab65332	100 test	Hemin Assay Kit	辅酶和辅助因子	六折
ab102504	100 test	Coenzyme A Assay Kit	辅酶和辅助因子	六折
ab204710	100 test	Flavin Adenine Dinucleotide (FAD) Assay Kit	辅酶和辅助因子	六折
ab204719	100 test	Acetate Assay Kit (Colorimetric)	辅酶和辅助因子	六折